

PDQuest™ 2-D Analysis Software

Version 6.2 Update Training Package

PDQuest software functionality and tools have been extensively enhanced. Additionally, a new focus on process automation and integration of data from a variety of sources will appeal to the proteome-oriented researcher. Highlights include:

- Greatly enhanced annotation capabilities that allow researchers greater flexibility and speed in annotating databases, as well as allowing spot annotation with URL addresses or document links to server path and file names
- More powerful data export and import capabilities
- More powerful image viewing capabilities and options, including multichannel viewing tools and multi-image reviewing tools
- Volume tools that allow researchers to override Gaussian spot quantitation
- Ability to build "cyber gels" with ReadyStrip™ narrow-range IPG strips
- Full integration with the ProteomeWorks™ system, including control of the Spot-Cutter excision robot as well as sample tracking through data analysis (with the Micromass™ MassPREP™ digestion robot and Micromass M@LDI™ and Q-TOF™ mass spectrometers)

Summary of major changes incorporated into PDQuest, version 6.2

Features/ Interface Design:

- Advanced Annotation Tool
- Improved Spot Data Export/Import
- Image Stack Tool
- Multichannel Image Merging
- Build Cyber Gels from Multiple Gels that Resolve Different pI or Molecular Weight Ranges
- Integrated ProteomeWorks Spot-Cutter Control
- Micromass Mass Spectrometry and Sample Prep Robot Integration
- The Scatter Plot Tool is linked to images, to allow identification of variant proteins
- Volume freehand and auto-contour tools have been added, allowing overwrite of Gaussian quantitation of spots
- Context-sensitive right-mouse-click menus
- Match-offset vectors now indicate outliers in red for better visibility when editing matching of spots

See subsequent pages of this training update for additional detail on the above features

Instrument Integration:

Bio-Rad Devices-

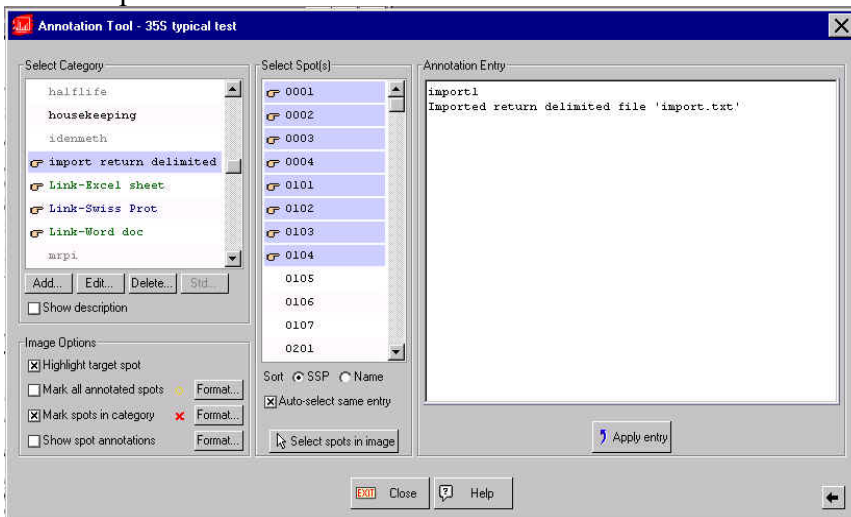
GSD-800 Scanning Densitometer
ProteomeWorks Spot Cutter

Other Devices-

Micromass mass spec data integration (MALDI & Q-TOF)

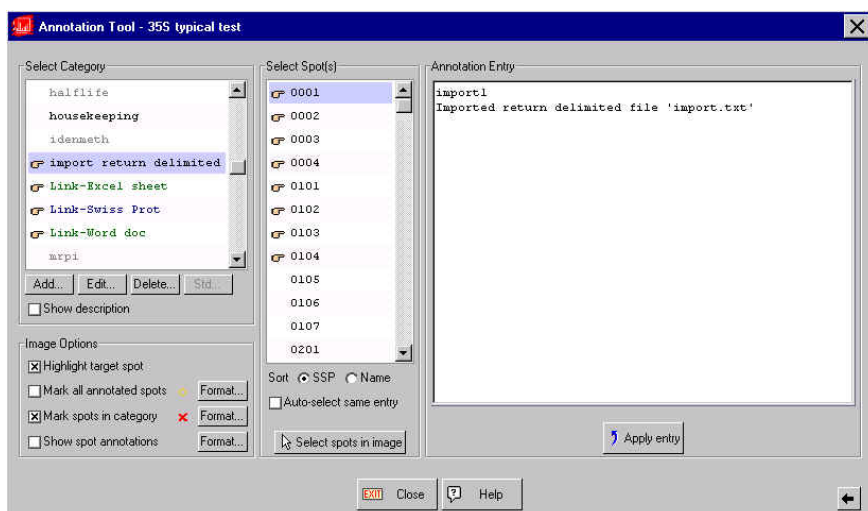
Advanced Annotation Tool

This tool is easier to work with; it stays resident until closed, and gives feedback on which categories have been annotated on a gel, and which spots within a category have been annotated. Multiple spots to be annotated can be selected from a gel or from an integral spot list and annotated in one operation. Conversion of entire analysis sets into annotations is now possible. You can also display multiple annotation categories simultaneously with unique, user-definable annotation markings. The Annotation Tool allows file path links to documents, including URL links, as annotations, with automatic launching of a web browser to the specified URL. It also dynamically builds HTML pages to view spot annotations.



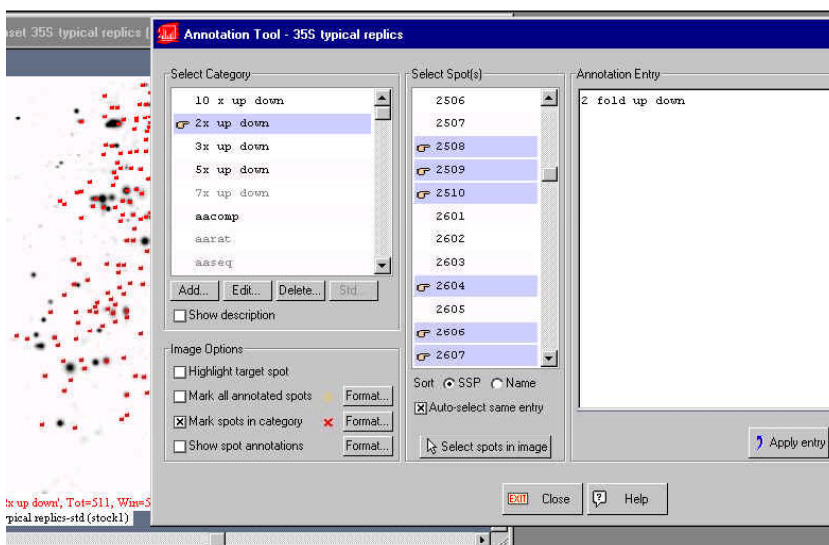
annotation 1.

- Positive indication of which categories have annotation entries.
- Positive indication of which spots within a category have annotation entries.
- Color-keyed indication of category format;
 - black for text annotation
 - blue for URL link annotation
 - green for file path name annotation (documents and spreadsheets)
- Positive indication of currently selected category and spot(s)
- Large field allows high annotation entry content.
- Enabling “Auto-select same entry” option allows all spots within a category with the same annotation entry to be selected and/or edited simultaneously in a single operation.



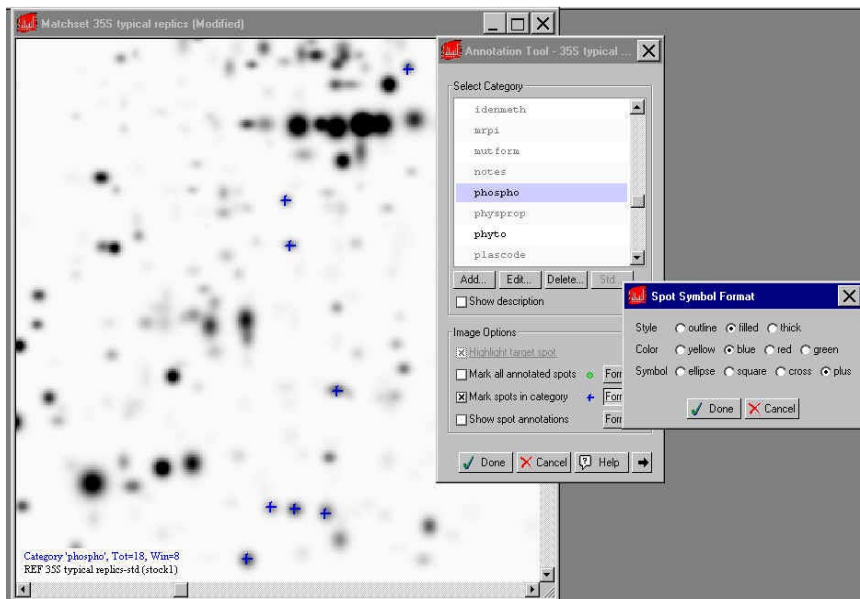
annotation 2.

- Disabling “Auto-select same entry” option allows selection or editing of a single spot within a category even if spots with same entry exist



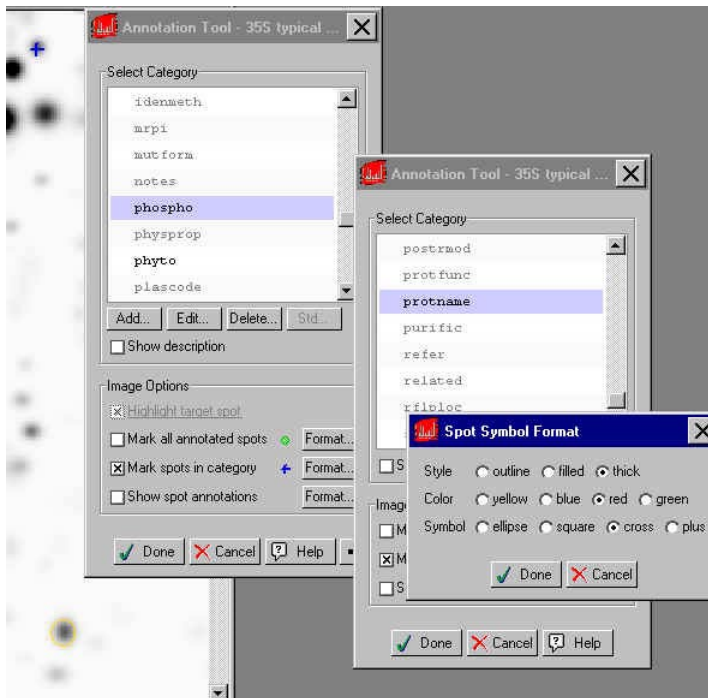
annotation 3.

- Can use Shift-click and Ctrl-click conventions to select multiple spots from spot list
- Can use Ctrl-click convention to select multiple spots from gel image



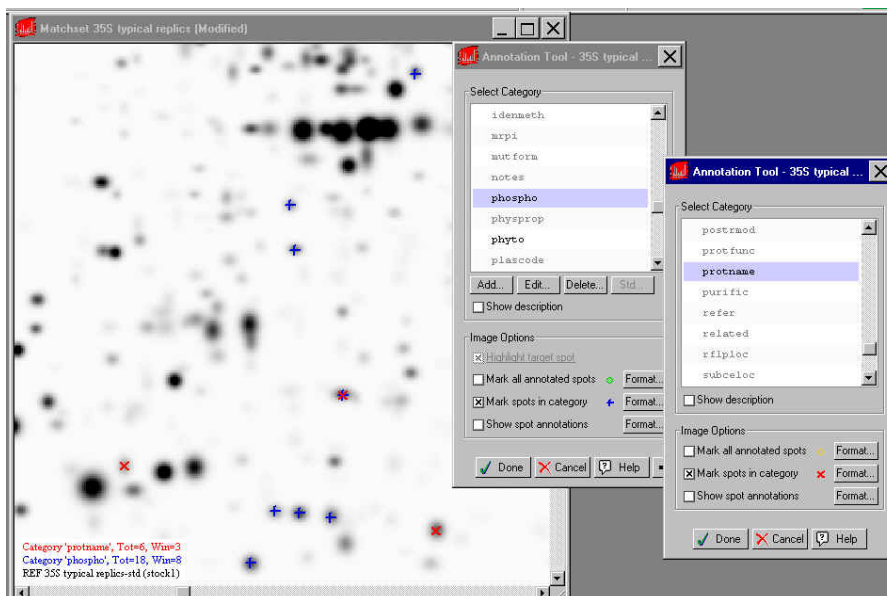
annotation 4.

- Gel image symbol overlay display options include;
 - Spots currently targeted (selected) for annotation entry
 - All annotated spots
 - Spots in selected category (*shown in above example*)
 - Annotation Category name
- User can define format of annotation overlay symbol for any specific category
 - style
 - color
 - symbol shape



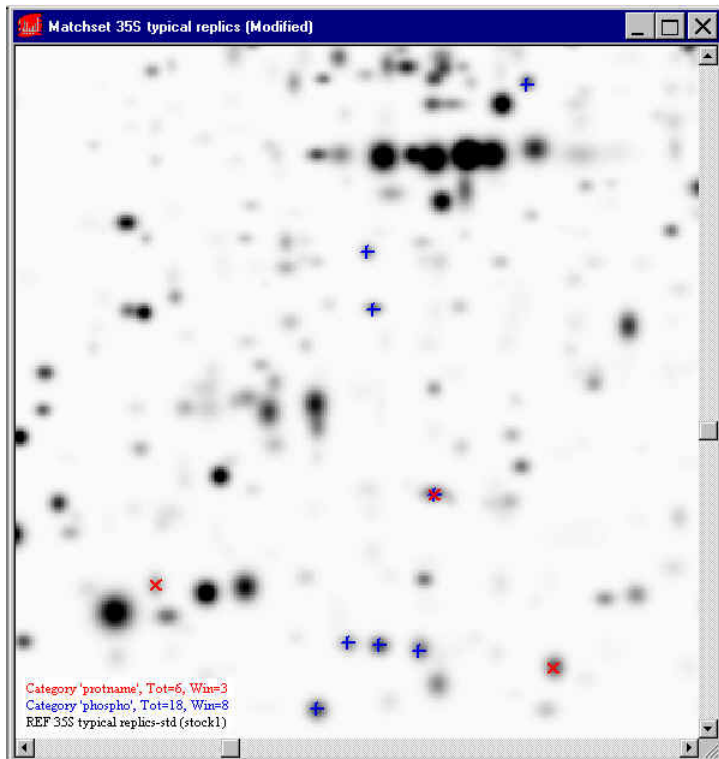
annotation 5.

- Tool stays resident until dismissed
- Multiple iterations of tool can be enabled, allowing simultaneous view of multiple categories
 - annotation overlay symbol can be defined independently for multiple categories



annotation 6.

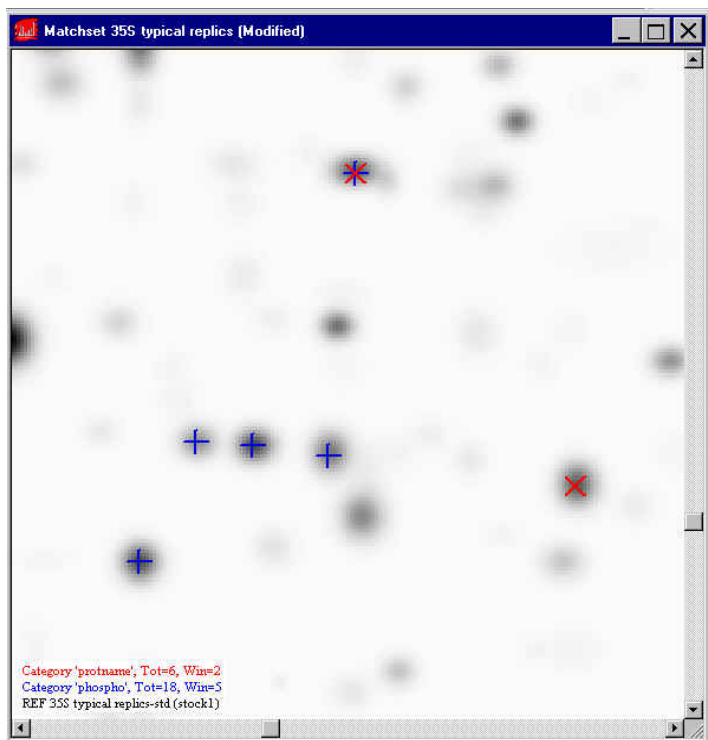
- Simultaneous view of two categories with unique spot symbol overlays
 - phosphorylase proteins (category phospho) as blue pluses
 - proteins with defined protein names (category protname) as red crosses



annotation 7.

CLOSE-UP VIEW

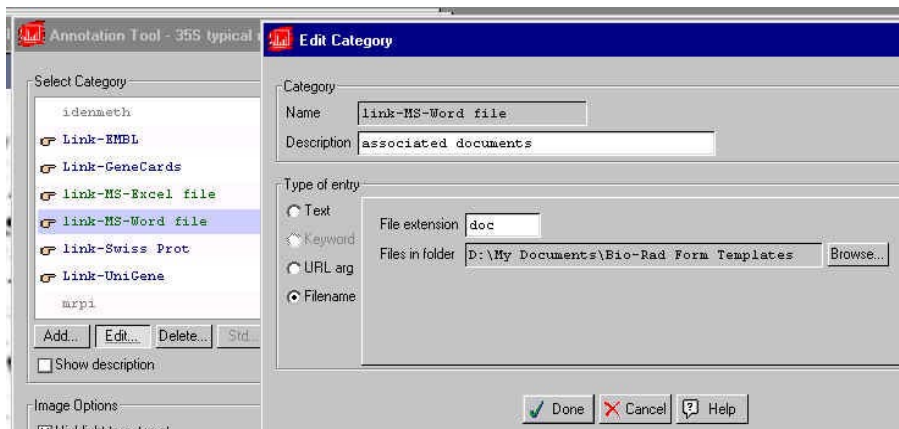
- Simultaneous view of two categories with unique spot symbol overlays
 - phosphorylase proteins (category phospho) as blue pluses
 - proteins with defined protein names (category protname) as red crosses
- Note category legend and spot counts in bottom left corner of window



annotation 8.

REALLY CLOSE-UP VIEW

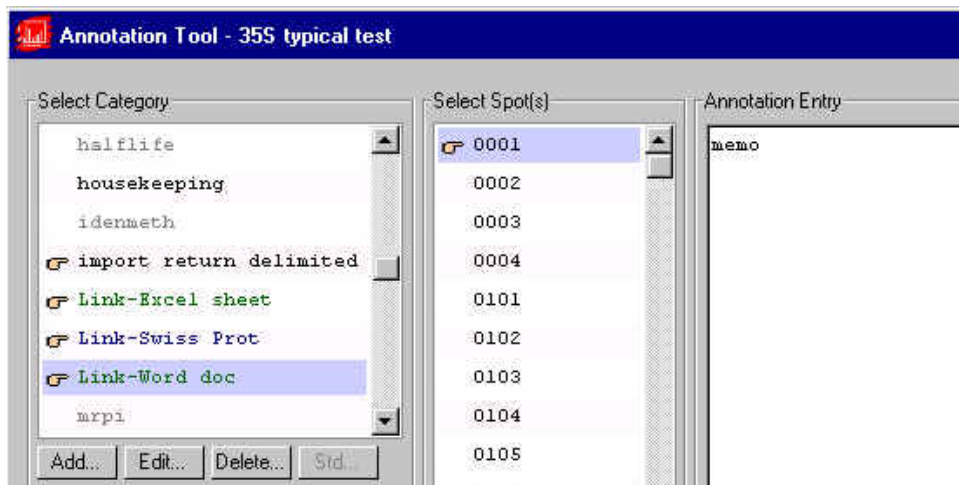
- Simultaneous view of two categories with unique spot symbol overlays
 - phosphorylase proteins (category phospho) as blue pluses
 - proteins with defined protein names (category protname) as red crosses
- Note category legend and spot counts in bottom left corner of window
- Note dual-symbol overlay for spot near top of window



annotation 9.

FORMATTING ANNOTATIONS: File Path Names

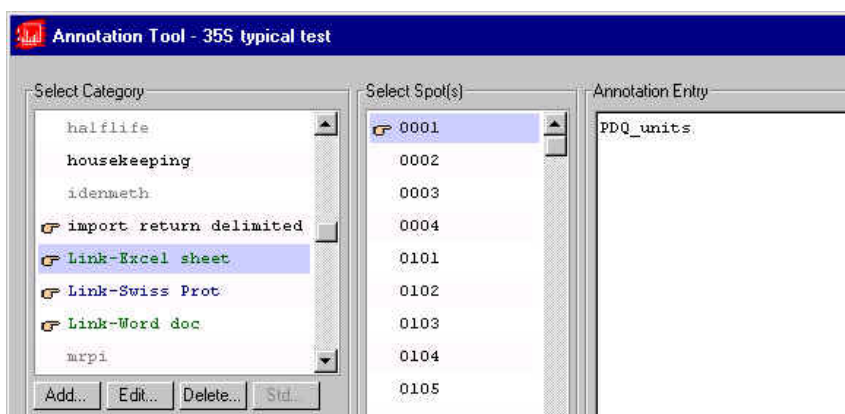
- File path links to MS-Word document or MS-Excel documents
- Automatic launch of application to link from right mouse menu
- File path and file name extensions are pre-defined for category



annotation 10.

ANNOTATION ENTRIES: File Path Names

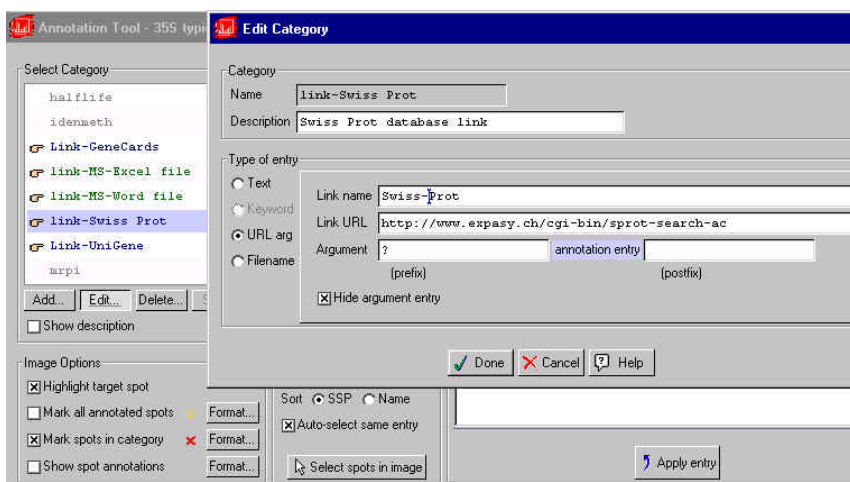
- File path link to MS-Word document
- File name extension not required



annotation 11.

ANNOTATION ENTRIES: File Path Names

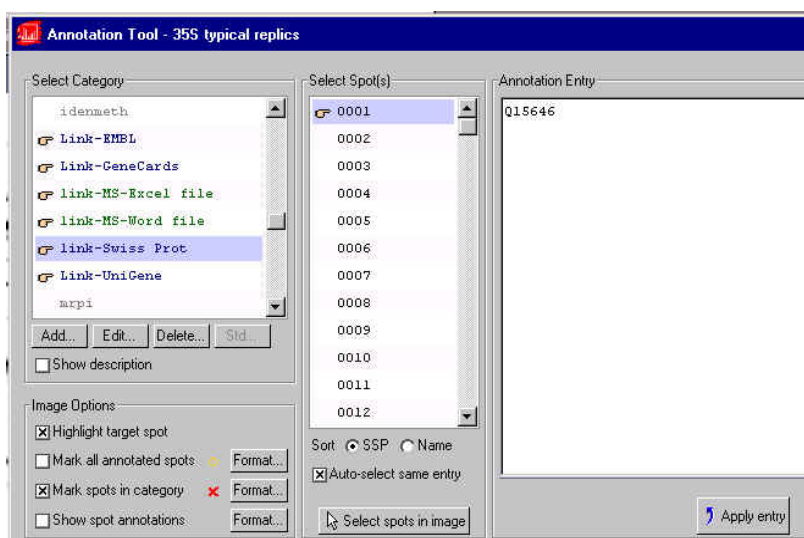
- File path link to MS-Excel document
- File name extension not required



annotation 12.

FORMATTING ANNOTATIONS: URL Links

- Link to URL
- Define URL link address
- Define argument prefix (to annotation entry)
- Define argument postfix (to annotation entry)



annotation 13.

ANNOTATION ENTRIES: URL Links

- Link to URL
- Annotation entry is database entry accession number

Spot information: SSP 1

unknown; Mr = 27.3 kDa ; pI = -1.0

Category	Description	Annotation Entry	Link
Link-EMBL	EMBL Database Link		EMBL
Link-GeneCards	Weizmann Inst GeneCards link		GeneCards
link-Swiss Prot	Swiss Prot database link		Swiss-Prot
Link-UniGene	NCBI UniGene database link		UniGene
10 x up down			
2x up down			
3x up down			
5x up down			
7x up down			
aacomp	Amino Acid Composition	AARTMR	
aarat	Amino Acid Ratios		

annotation 14.

BROWSE ANNOTATIONS PAGE:

- HTML page view of spot annotation
- Launched from right-mouse menu from an annotated spot
- Dynamically built in local browser to show all spot annotation updates
- Shows all categories and presence of annotation entries
- Shows File Path Name links
- Shows URL links as hyperlinks

ExPASy Home page	Site Map	Search ExPASy	Contact us	SWISS-PROT
Mirror sites: Australia Canada China Taiwan				

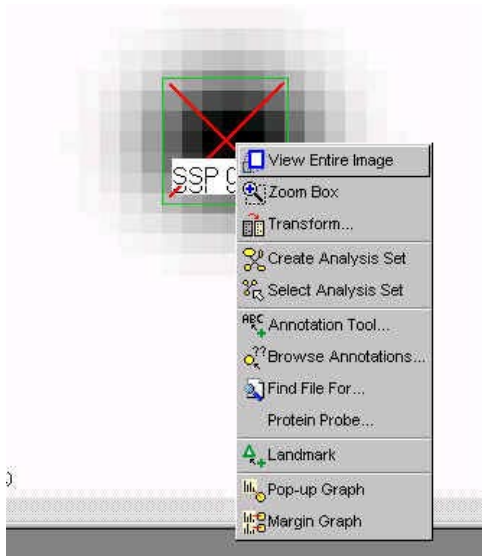
NiceProt View of SWISS-PROT: [Q15646](#)

[\[General\]](#) [\[Name and origin\]](#) [\[References\]](#) [\[Comments\]](#) [\[Cross-references\]](#) [\[Keywords\]](#) [\[Features\]](#) [\[Sequence\]](#) [\[Tools\]](#)

General information about the entry	
Entry name	25AL_HUMAN
Primary accession number	Q15646
Secondary accession number(s)	O75686
Entered in SWISS-PROT in	Release 35, November 1997
Sequence was last modified in	Release 39, May 2000
Annotations were last modified in	Release 39, May 2000
Name and origin of the protein	
Protein name	59 KDA 2'-5' OLIGOADENYLATE SYNTHETASE LIKE PROTEIN
Synonym(s)	EC 2.7.7.- P59OASL THYROID RECEPTOR INTERACTING PROTEIN 14 TRIP14
Gene name(s)	OASL OR TRIP14
From	Homo sapiens (Human)
Taxonomy	Eukaryota; Metazoa; Chordata; Craniata; Vertebrata; Euteleostomi; Mammalia; Eutheria; Primates; Catarrhini; Hominidae

annotation 15.

- Link to URL destination from BROWSE ANNOTATIONS PAGE:
- Automatic launch of local Web Browser



annotation 16.

- Open linked documents through right-mouse menu from spot
 - Use “Find File For...” command
 - Automatic launch of local application
-

Improved Spot Data Export/Import

Export of all matchset data is now possible. You can import a list of SSP numbers as a file, and convert it into an analysis set.

Export Matchset - Matchset 355 typical replics

Data to export

Spot data by gel | Spot data by group

☒ Center position ☒ Quantity ☒ Quality

☒ Gauss size ☒ Norm. Quantity ☒ Peak value

Saturated spots: ☒ Estimate ☐ (-3.0) ☐ Other:

Unmatched spots: ☒ Estimate ☐ (-1.0) ☐ Other:

☒ SSP ☐ DSN ☐ Group Id ☒ Mr. pl

Annotations: ☐ None ☐ All ☒ Specified:

Analysis sets: ☐ None ☒ All ☐ Specified

Spots to export

☐ All spots

☐ Matched to all gels

☒ In analysis set:

Spot count:

Output options

☒ Write column headers

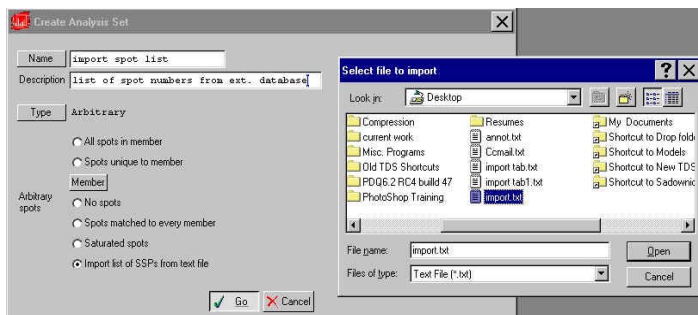
Sort by: ☒ SSP ☐ DSN ☐ Gel/group

Separators: ☒ Tabs [Excel] ☐ Commas [1-2-3]

Export to: ☐ File ☒ Clipboard

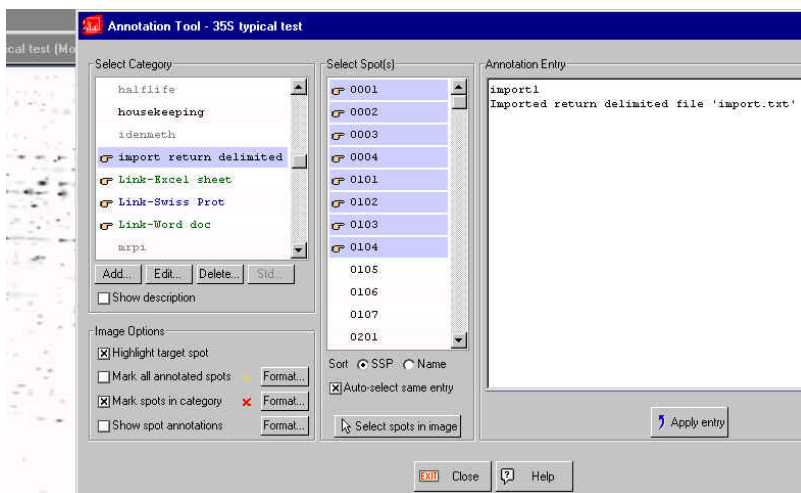
export 1.

- Check boxes allow tailoring of any or all parameters to be included in export file
- Parameters can include annotation category entries and/or analysis set membership status
 - options include none, all , or specified category or analysis set
- Can restrict selection of spots for export;
 - all spots
 - spots matched to all gels
 - spots in a specific analysis set
- Export to clipboard allows convenient paste into spreadsheet programs for immediate viewing



import 1.

- Create an Arbitrary Analysis Set
- Define Type as “Import List of SSPs from text file”
- Browse to return-delimited text file containing spot list
-valid spot entries are “101” or “0101”

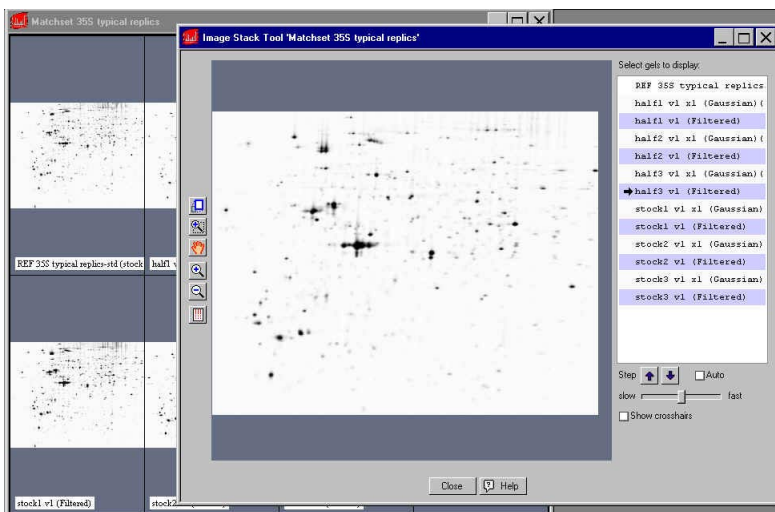


import 2.

- Convert Analysis Set into an Annotation Category
-from menu: **Analysis > More Annot. Tools > Create Annotation from Set**

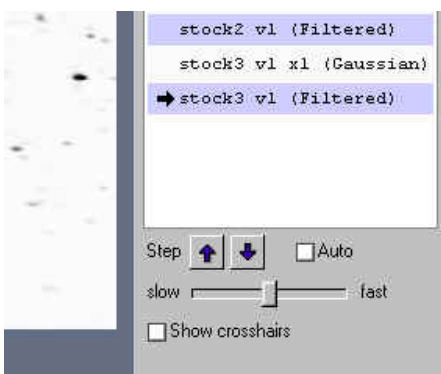
Image Stack Tool

This tool allows aligned-stack viewing of multiple gels, and also offers an auto-flicker option with user-adjustable speed. Aligned stack viewing is only available if gels have been matched. Local alignment will be performed within any window view based on the spot(s) closest to the center of the view.



stack 1.

- All open images are loaded into image list window
- Use Shift-Click and Ctrl-click conventions to choose images for viewing
- If matched

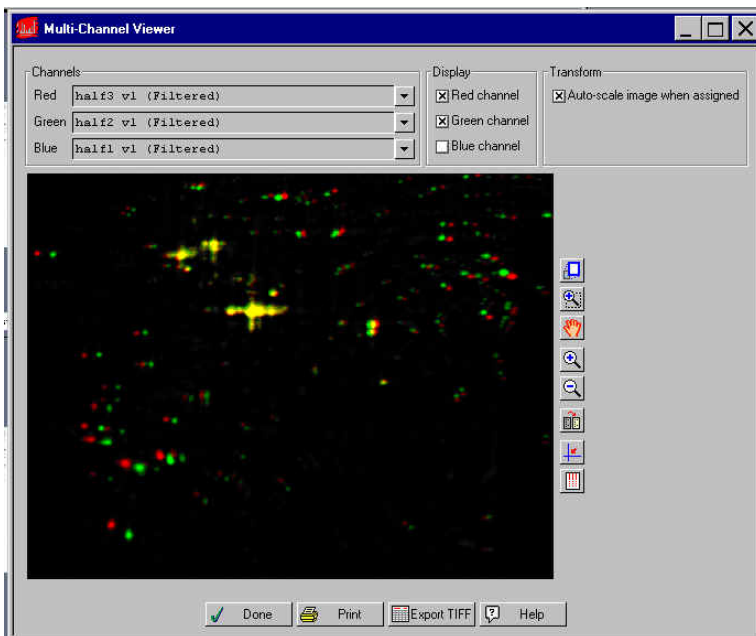


stack 2.

- flicker tools
 - up/down
- auto-flicker tools
 - on/off
 - speed of image flickering

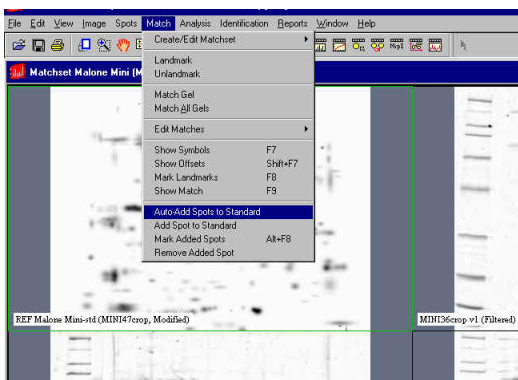
Multichannel Image Merging

Merge up to 3 image channels with independent red, green, or blue color display. Balance the display with independent channel transformation. Zoom in to see fine details like spot overlap.



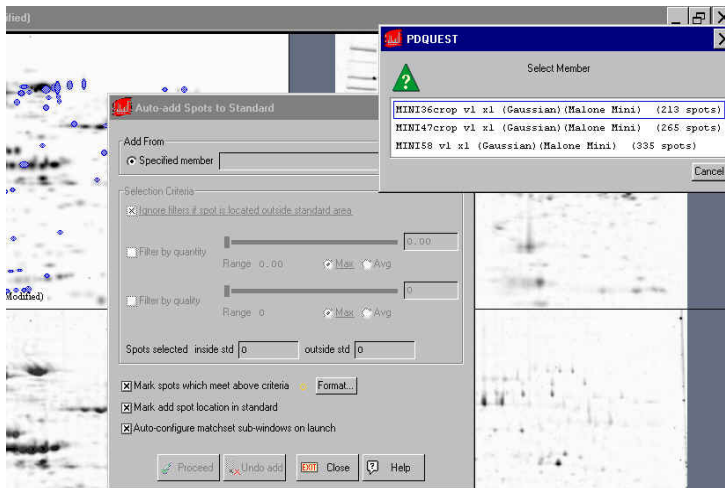
Build Cyber Gels from Multiple Gels that Resolve Different pI or Molecular Weight Ranges

The new Auto-add to Standard function allows addition of all unique spots from a given gel to the standard reference image in one operation. These tools allow expansion of the standard reference image to accommodate a larger area (for example, when combining 2 different IPG strip separations) to build a cyber gel.



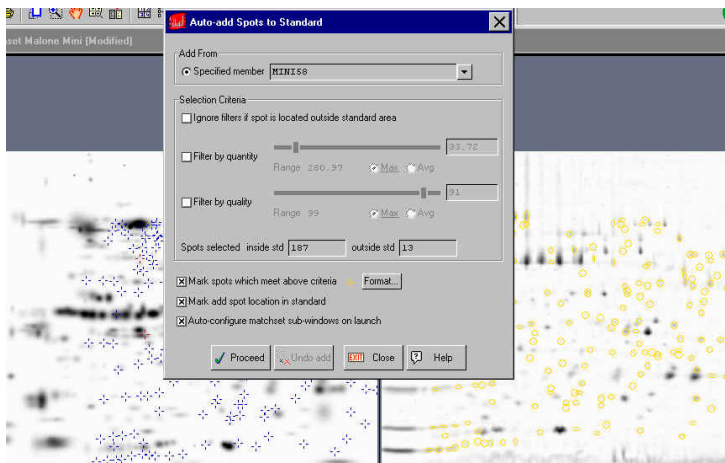
Cyber 1.

- Menu : **Match > Auto-add to standard**



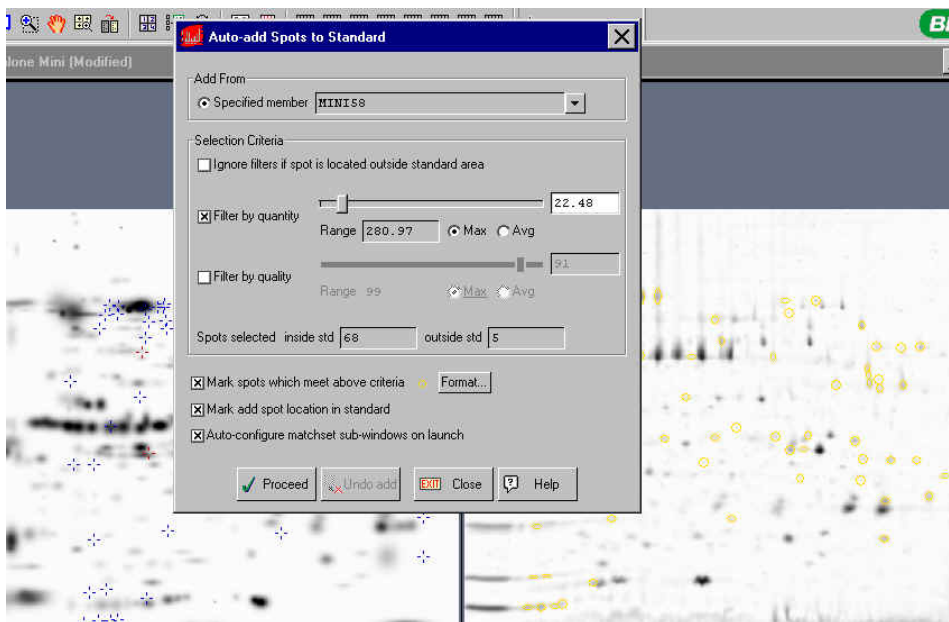
Cyber 2.

- Choose gel to add unmatched spots from



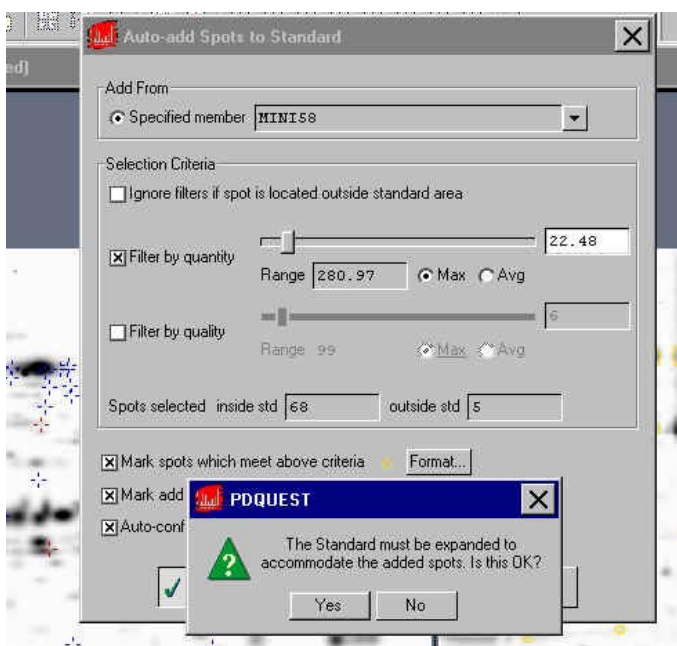
Cyber 3.

- Gel image overlay with targeted spots in member and targeted position to add in Standard Reference



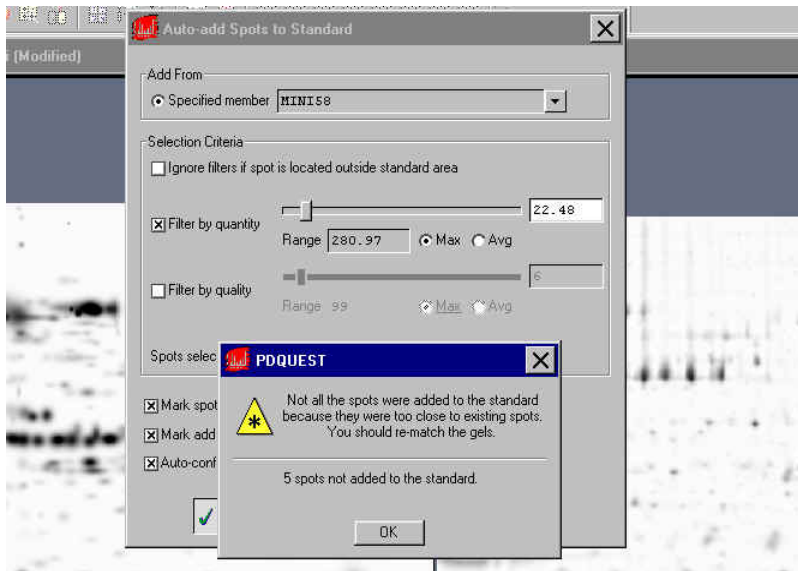
Cyber 4.

- Filtering of spot selection by both **quantity** and **quality**
- Instant feedback in gel image and spot selection number



Cyber 5.

- Ability to expand Standard Reference to accommodate a wider pH or MW range
- User can select to ignore filters if targeted spot is outside the Standard Reference area

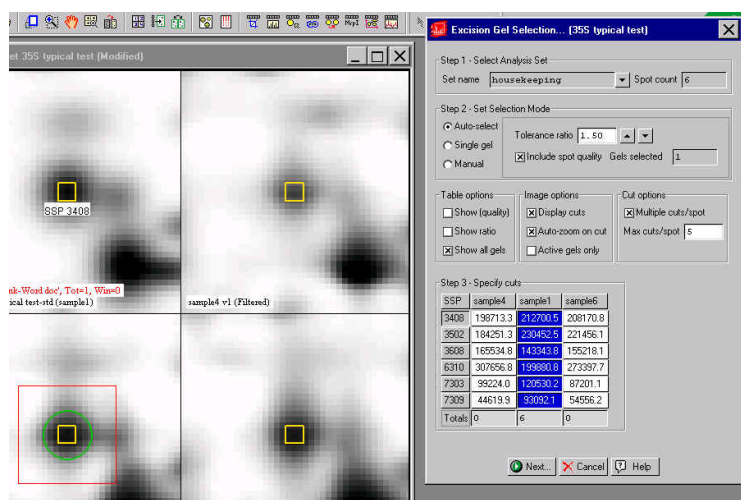


Cyber 6.

- Proximity over-ride and warning

Integrated ProteomeWorks Spot-Cutter Control

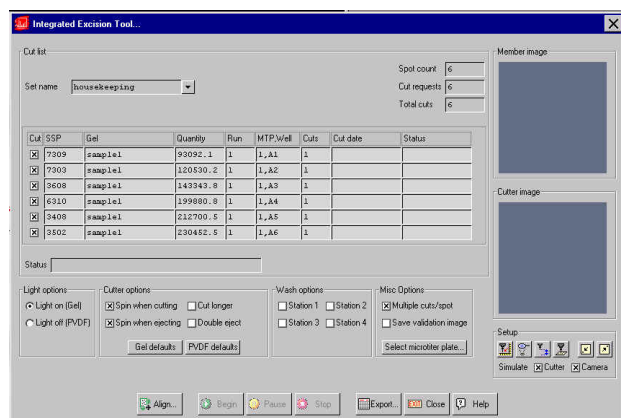
The software will drive the spot cutter to cut out any analysis set defined in PDQuest. In PDQuest software, with automatic targeting of spots to the correct location on the gel. A Basic Excision (point-and-click) cutting mode is also available. Basic Excision will also run in PDQuest Viewer mode.



Spot Cutter 1.

GEL SELECTION TOOL

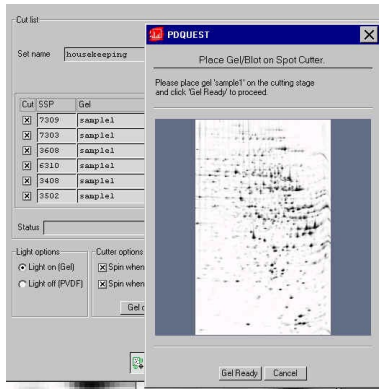
- Spots are pre-selected from PDQuest Analysis Sets
- Spot Cutter software can automatically target the gel with the highest expression of any given spot (to maximize excision yield)
 - automatically tracks and prompts for correct gel during excision
 - user can over-ride or supplement automatic selection
- Targeted spots in gel images presented for easy verification



Spot Cutter 2.

INTEGRATED EXCISION TOOL

- Controls alignment and cutting/re-cutting operations



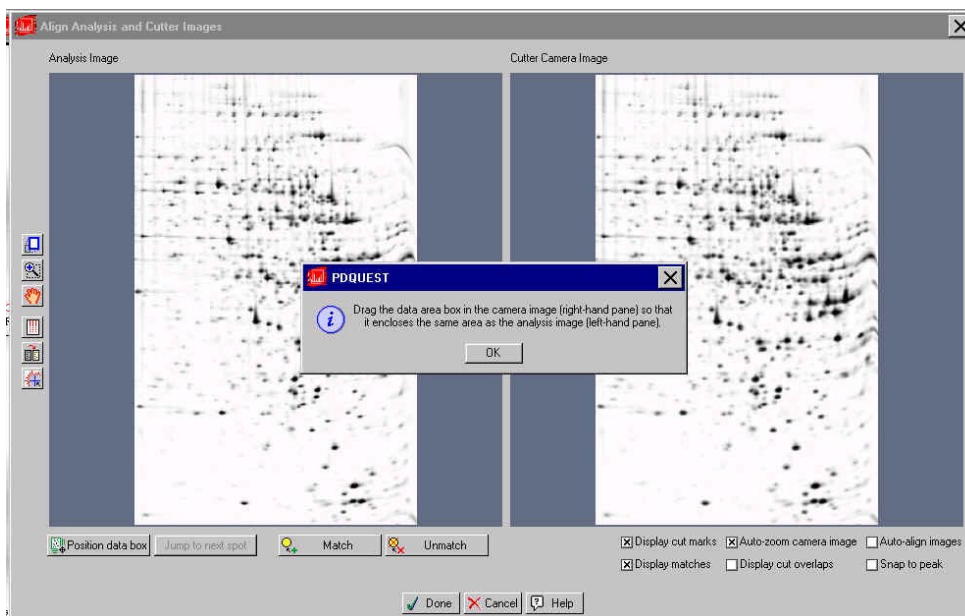
Spot Cutter 3.

- Prompt for appropriate gel



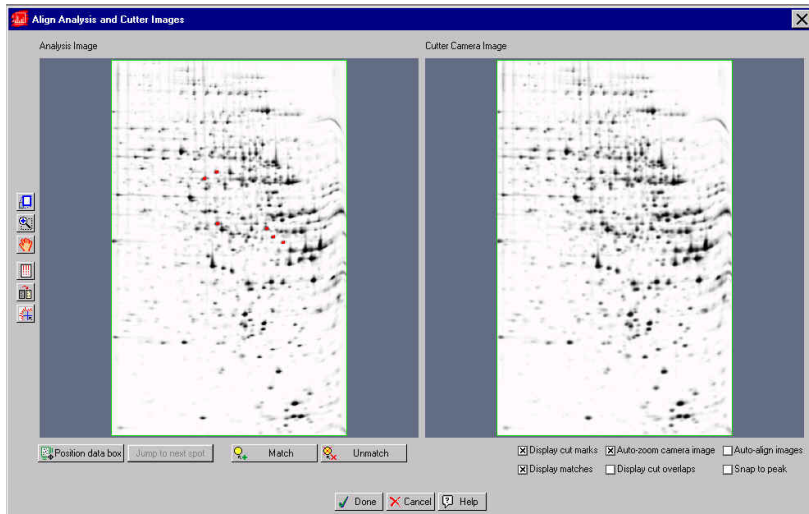
Spot Cutter 4.

- Spot Cutter Simulation mode for demos/training



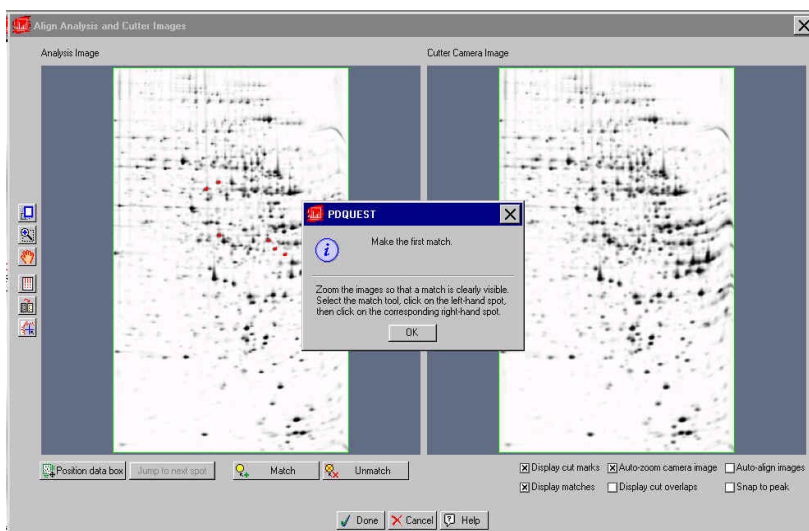
Spot Cutter 5.

- Prompts for initial alignment; gel area



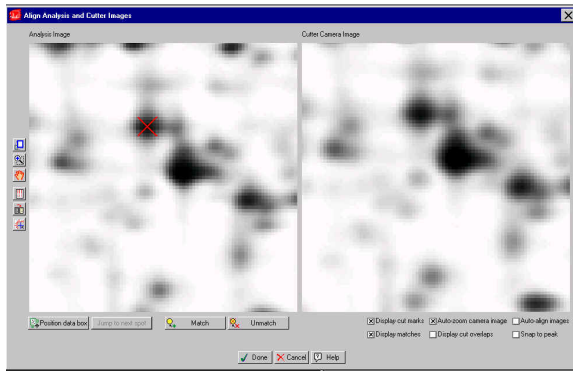
Spot Cutter 6.

- gel area aligned



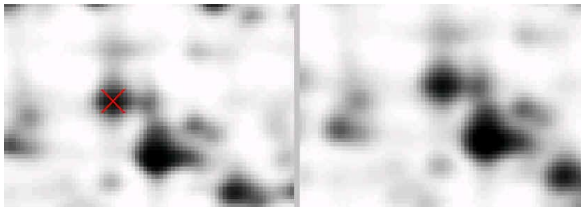
Spot Cutter 7.

- Prompts for gel alignment; targeted spot in analysed image versus gel on Spot Cutter table



Spot Cutter 8.

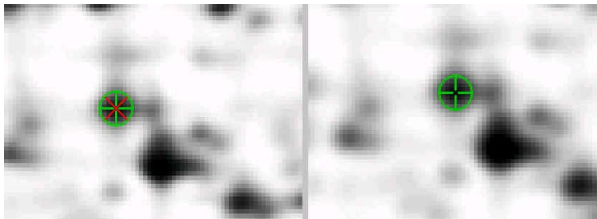
- First targeted cut
 - Analysis image on left, Spot Cutter target on image on right
-



Spot Cutter 9.

CLOSE UP

- First targeted cut
 - User clicks on matching spot in right side image to confirm cut location
-



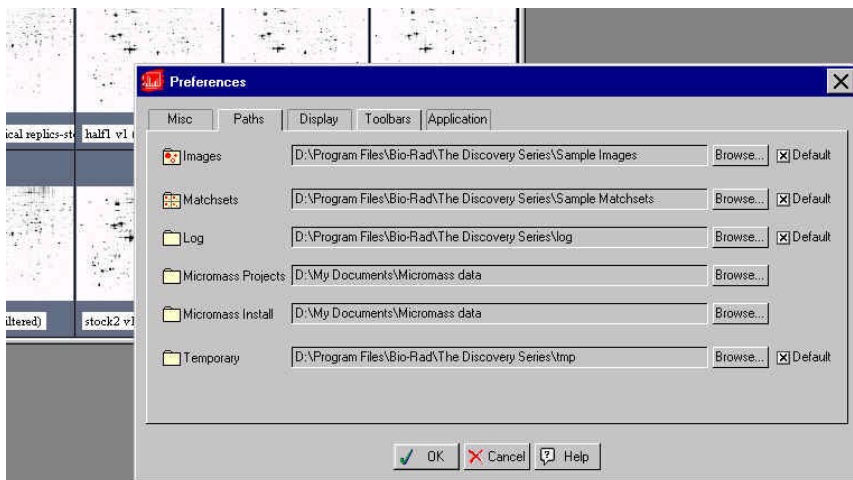
Spot Cutter 10.

CLOSE UP

- First targeted cut after alignment
- Target confirmation can be dragged to adjust accuracy of cutter tip placement
- Green circle is actual size of cutting tip
- After alignment, Shift-Click to jump to next targeted cut

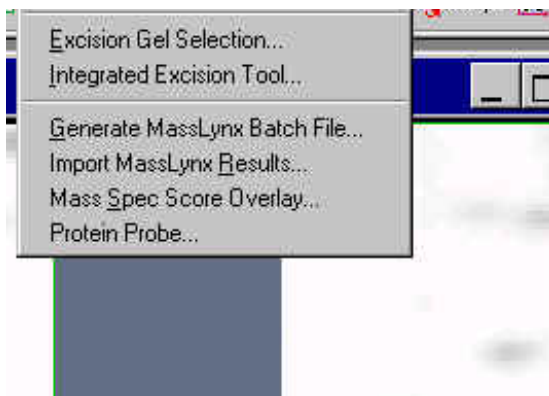
Micromass Mass Spectrometry and Sample Prep Robot Integration

You can track the spot (after excision with the spot cutter) through the trypsin digestion robot and to the data file generated for that spot. Automatically update all spot annotations in a matchset with protein identification annotations from the resulting Micromass data files. Launch Micromass software via the ProteinProbe(tm) browser to see the associated mass spectra for a spot.



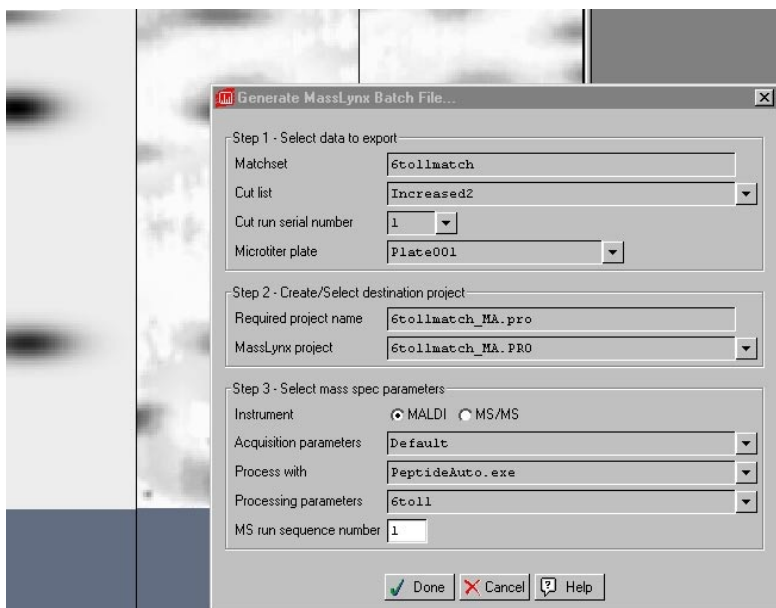
Micromass 1.

- Define pathways for Micromass data exchange
 - where to export data file to establish a MassLynx project file
 - where to find resulting protein identification results file



Micromass 2.

- Export a batch file which contains the necessary info to establish links to the MassLynx project name for the mass spec results and instrument / processing software control parameters
- Import the resulting data file with automatic annotation category creation and data entry for identified spots



Micromass 3.

- Export file defines MassLynx project name and sets Mass Spec parameters (for either MALDI or Q-TOF instruments)
 - Mass spec acquisition parameters
 - spectra processing parameters
 - database search/comparison parameters
- MassLynx project name must match PDQuest required name and must use default file name extensions

MS	Sample	File Name	File Text	MS File	Process	Parameter File
1	A.1	Increased2_1_Plate001_1_4702	Default	PeptideAuto.exe	bw01	
2	A.2	Increased2_1_Plate001_1_5101	Default	PeptideAuto.exe	bw01	
3	A.3	Increased2_1_Plate001_1_6801	Default	PeptideAuto.exe	bw01	
4	A.4	Increased2_1_Plate001_1_4801	Default	PeptideAuto.exe	bw01	
5	A.5	Increased2_1_Plate001_1_3901	Default	PeptideAuto.exe	bw01	
6	A.6	Increased2_1_Plate001_1_7401	Default	PeptideAuto.exe	bw01	
7	A.7	Increased2_1_Plate001_1_2201	Default	PeptideAuto.exe	bw01	
8	A.8	Increased2_1_Plate001_1_0301	Default	PeptideAuto.exe	bw01	
9	A.9	Increased2_1_Plate001_1_4603	Default	PeptideAuto.exe	bw01	
10	A.10	Increased2_1_Plate001_1_0202	Default	PeptideAuto.exe	bw01	
11	A.11	Increased2_1_Plate001_1_5301	Default	PeptideAuto.exe	bw01	
12	A.12	Increased2_1_Plate001_1_7302	Default	PeptideAuto.exe	bw01	
13	B.1	Increased2_1_Plate001_1_6302	Default	PeptideAuto.exe	bw01	
14	B.2	Increased2_1_Plate001_1_0801	Default	PeptideAuto.exe	bw01	
15	B.3	Increased2_1_Plate001_1_2202	Default	PeptideAuto.exe	bw01	
16	B.4	Increased2_1_Plate001_1_6804	Default	PeptideAuto.exe	bw01	
17	B.5	Increased2_1_Plate001_1_5402	Default	PeptideAuto.exe	bw01	
18	B.6	Increased2_1_Plate001_1_7701	Default	PeptideAuto.exe	bw01	
19	B.7	Increased2_1_Plate001_1_7801	Default	PeptideAuto.exe	bw01	
20	B.8	Increased2_1_Plate001_1_2906	Default	PeptideAuto.exe	bw01	
21	B.9	Increased2_1_Plate001_1_1701	Default	PeptideAuto.exe	bw01	
22	B.10	Increased2_1_Plate001_1_4701	Default	PeptideAuto.exe	bw01	
23	B.11	Increased2_1_Plate001_1_6701	Default	PeptideAuto.exe	bw01	
24	B.12	Increased2_1_Plate001_1_7901	Default	PeptideAuto.exe	bw01	
25	C.1	Increased2_1_Plate001_1_2801	Default	PeptideAuto.exe	bw01	
26	C.2	Increased2_1_Plate001_1_3303	Default	PeptideAuto.exe	bw01	
27	C.3	Increased2_1_Plate001_1_7902	Default	PeptideAuto.exe	bw01	
28	C.4	Increased2_1_Plate001_1_4902	Default	PeptideAuto.exe	bw01	

Micromass 4.

- Project file is set up within MassLynx software by mass spec operator

Import MassLynx Results...

Step 1 - Select project

Matchset: 6tollannotated2

Instrument: ☒ MALDI ☐ MS/MS

Required project name: 6tollannotated2_MA.pro

MassLynx project: 6tollmatch_MA.PRO

Step 2 - Select run

Step 3 - Select sample

Sample	Date
E,7	31-AUG-2000
E,8	31-AUG-2000
E,9	31-AUG-2000
E,10	31-AUG-2000
E,11	31-AUG-2000

Step 4 - Select primary hit

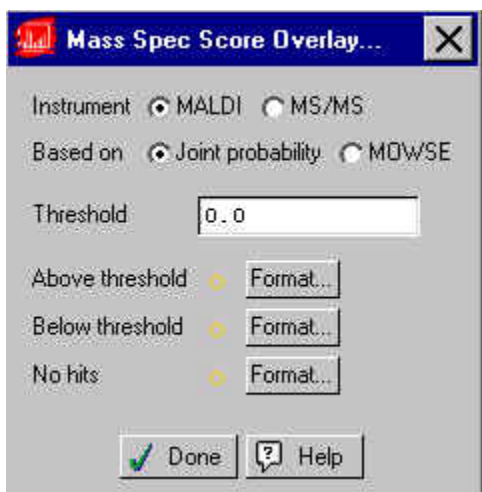
Gene	Prot	Peptides	Coverage	MW	pI	Score
[gene=tolB]	[prot=TolB protein]	21 peptides	55.79% coverage	47.72kDa	9.07pH	score 119657.9
[gene=PA3803]	[prot=conserved hypothetical protein]	6 peptides	18.87% coverage	40.06kDa	6.38pH	score 6579.1
[gene=pncB1]	[prot=nicotinate phosphoribosyltransferase]	5 peptides	14.79% coverage	46.10kDa	6.26pH	score 4904.2
[gene=PA3067]	[prot=probable transcriptional regulator]	4 peptides	27.21% coverage	16.37kDa	10.07pH	score 4883.3
[gene=tig]	[prot=trigger factor]	5 peptides	14.22% coverage	48.58kDa	4.87pH	score 4297.1
[gene=trmA]	[prot=tRNA (uracil-5-)-methyltransferase]	4 peptides	16.80% coverage	41.21kDa	6.89pH	score 3781.3
[gene=PA3343]	[prot=hypothetical protein]	4 peptides	8.74% coverage	43.60kDa	8.29pH	score 3549.7

Score: ☒ Joint probability ☐ MOWSE

Buttons: Apply, Protein Probe..., EXIT, Close, Help

Micromass 5.

- File import after MS run complete includes “most probable hit selection” dialogue
- Hit data summary for probable hits includes
 - number of matched peptides in spectra
 - percentage of the targeted protein covered by the identified peptides
 - molecular weight and isoelectric point of the probable match
 - match probability score (user selectable for Joint Probability or MOWSE algorithms)



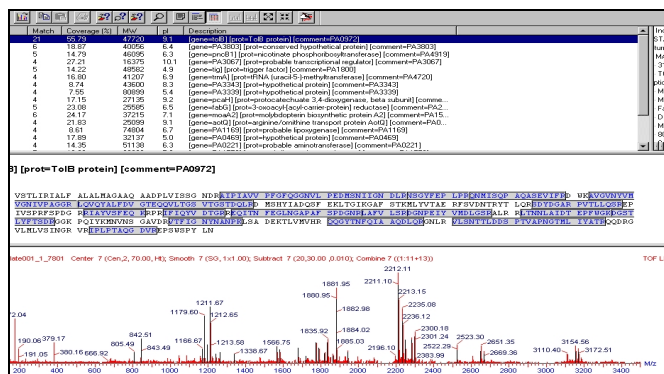
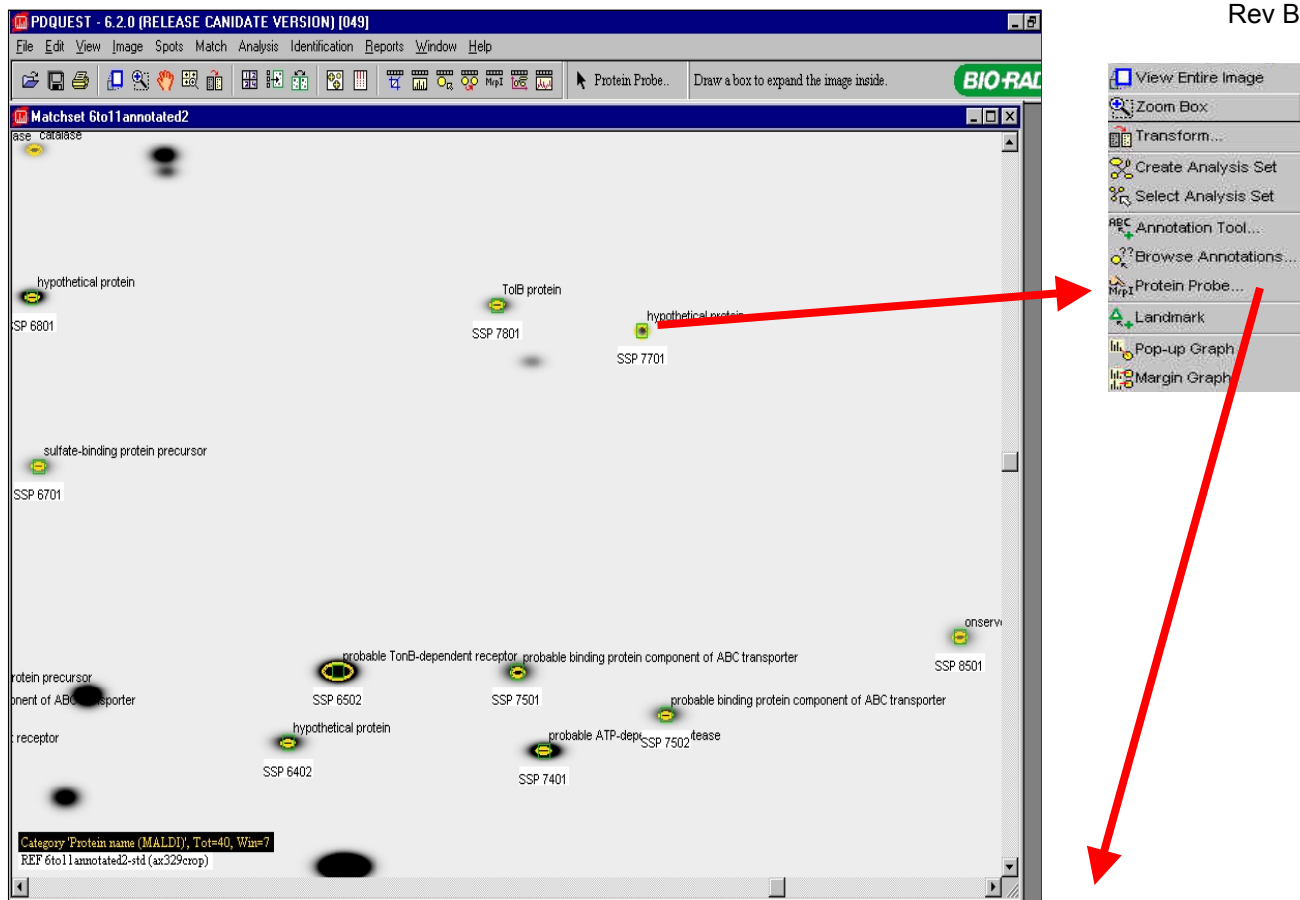
Micromass 6.

- Identified proteins in the gel image can be highlighted with hit indicator overlays based on the Match Probability Score
- Overlays can be color coded for user-defined confidence thresholding



Micromass 7.

- Import MassLynx data files to automatic create Annotation Categories and/or update Annotation Category entries with protein identification info

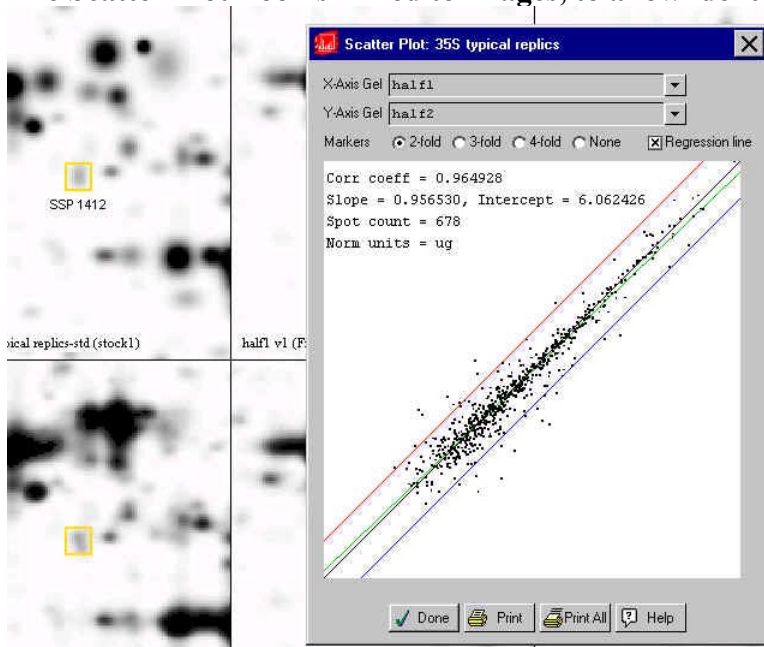


Micromass 8.

- Launch ProteinProbe software to view actual mass spectra and structure information associated with protein identification
- Advantages of integrated 2D Analysis software / Spot Cutter control / Mass Spec data
 - “Eliminates errors”
 - “Cuts time by 80%”
 - “Allows gel map to be used as point of reference for other data - visual image”
 - “Facilitates user compliance”
 - “Allows study of more complex proteomes”

*ProteomeWorks system evaluation by Dr. Brad Walsh, APAF
September 2000, Sienna Proteome Conference*

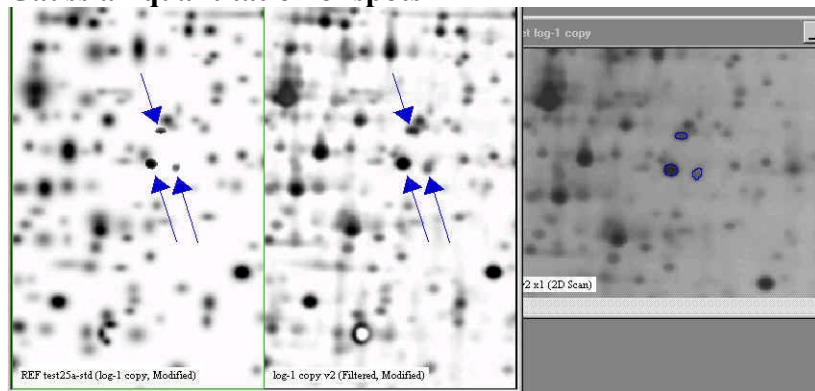
The Scatter Plot Tool is linked to images, to allow identification of variant proteins



scatter plot 1.

- Clicking on any data point in the plot will target the correlating protein on the gel
- Spots falling outside the standard deviation can now be easily identified on the gel image
- only matched spots are plotted in the tool

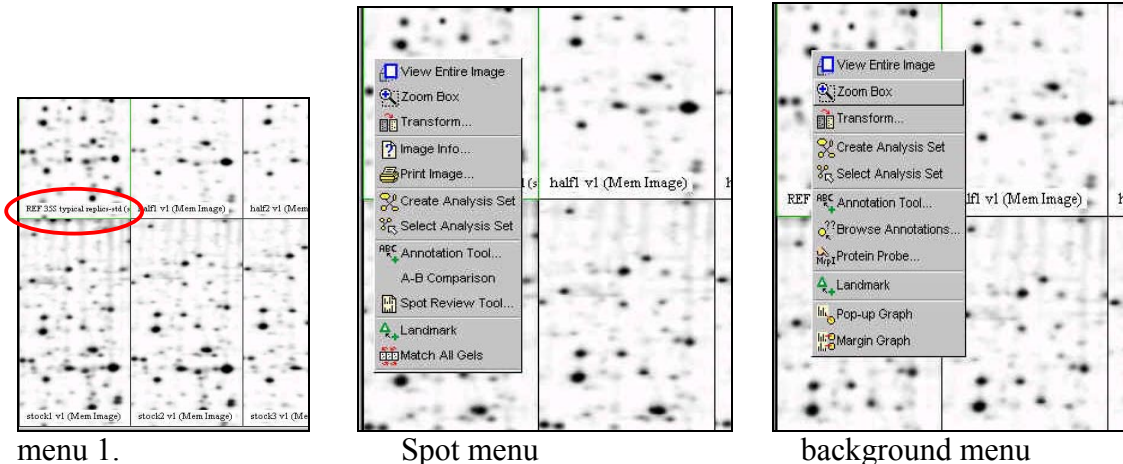
Volume freehand and auto-contour tools have been added, allowing overwrite of Gaussian quantitation of spots



contour 1.

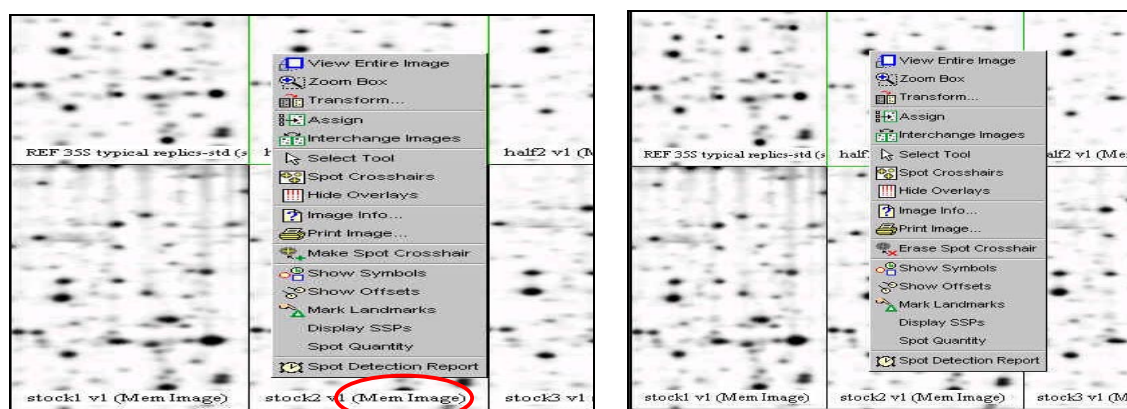
- Can apply a contour to scanset or matchset images

Context-sensitive right-mouse-click menus



Menus available from the **Standard Reference image**

-context referenced to cursor location on spots or gel background



Menus available from the **member images**

-context referenced to cursor location on spots or gel background

Match-offset vectors now indicate outliers in red for better visibility when editing matching of spots



offset 1.

- Offsets are compared to local neighbors for direction and length of distortion
- outliers are indicated in red

Ordering Information

Catalog # Description

- | | |
|----------|------------------------------------|
| 170-8603 | PDQuest 2-D Analysis Software, PC |
| 170-8611 | PDQuest 2-D Analysis Software, Mac |
| 170-8626 | PDQuest Software Upgrade, PC/Mac |