

ChemiDoc™ Touch Imaging System

Sensitivity in detection, power in quantitation



ChemiDoc Touch Imaging System

Best-in-class performance

Superior to film in signal-to-noise ratio

Equal to film in sensitivity and resolution

High-quality imaging of gels and western blots

Highly intuitive Image Lab[™] Touch Software

Streamlined path from experiment to usable data

Stain-free enabled

Publication-quality images at your fingertips

HIGH-PERFORMANCE IMAGING

EASY, FLEXIBLE INTERACTION

STAIN-FREE ENABLED

WESTERN BLOTTING CONSUMABLES

High-Performance Imaging Easy Acquisition Features As sensitive as film, with advanced blot detection technology to Includes image preview, auto-focus, auto-exposure, and additional determine best exposure for faint exposure options and intense bands **Convenience in Storing** and Sharing Data Export images via USB or network connection BIO RAD (6) Assess Images at the Point of Acquisition Smart Tray Technology™ Pinch and zoom images on the 12-inch Automatically recognizes touch screen; access a range of tools your application with Image Lab Touch Software GelGreen or any Chemiluminescent Coomassie Blue, silver, SYBR® Stains. blots, stain-free gels/ and other stains. blots, and ethidium bromide, SYPRO Ruby, and other stains.

HIGH-PERFORMANCE IMAGING

Get the sensitivity of film without the hassles of film processing, darkroom chemicals, or associated mishaps. Combine this sensitivity with a suite of tools to optimize imaging and quantitation, and achieve an unmatched ability to resolve the faintest and most intense bands into meaningful data.

The ChemiDoc Touch Imaging System is comparable to film ...

Detect low signal at the same exposure time

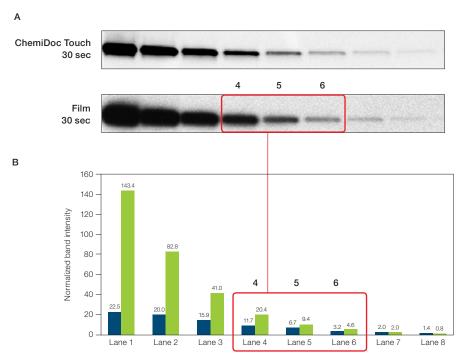


Fig. 1. Comparison of sensitivity between the ChemiDoc Touch Imaging System and film. A, western blot analysis of Lacl expression was conducted using 2x serial dilutions (starting at 0.31 µg protein) of E. coli cell lysate. The membranes were either imaged on the ChemiDoc Touch Imaging System for 30 sec or exposed to film for 30 sec. B, the normalized band densities illustrate the ability of the ChemiDoc Touch Imaging System to detect low signal bands at the same exposure time as film. The red boxes represent the limited linear dynamic range of film. ChemiDoc Touch Imaging System, 30 sec (■); film, 30 sec (■).

... and in many cases the ChemiDoc Touch Imaging System is superior to film.

Reveal faint protein bands missed by film.

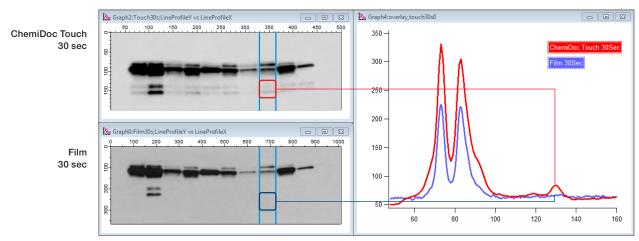


Fig. 2. Side-by-side comparison between the ChemiDoc Touch Imaging System and film. Levels of the three isoforms of the pro-apoptotic protein Bim were measured in various cell lines using western blot analysis. The membranes were either imaged on the ChemiDoc Touch Imaging System for 30 sec or exposed to film for 30 sec to compare detection sensitivities. As shown by the overlay graph, the ChemiDoc Touch Imaging System was better able to detect faint protein bands than film.



Best-in-Class Digital Image Quality

Comparison of the ChemiDoc Touch Imaging System with other digital imagers

A
ChemiDoc Touch Imaging System — 15 sec exposure



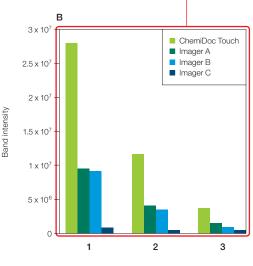
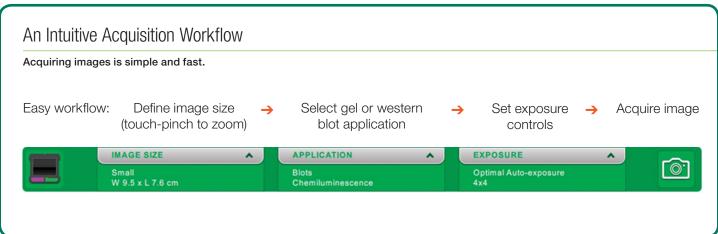
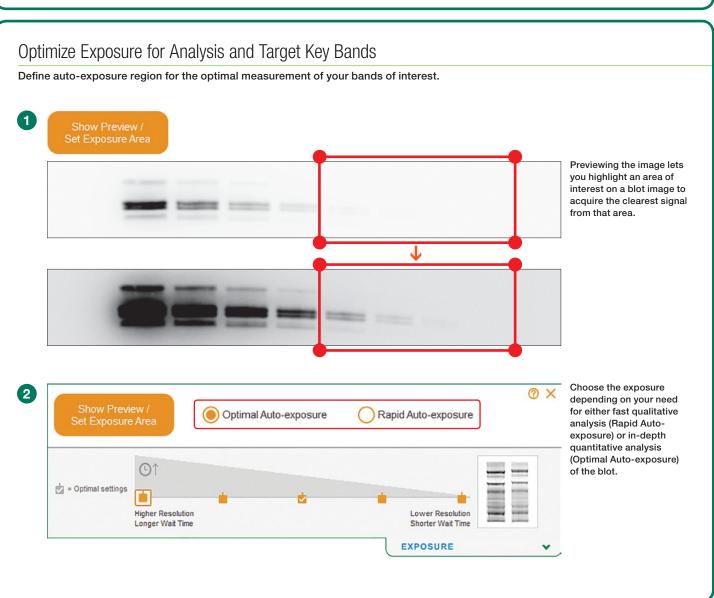


Fig. 3. Comparison between the ChemiDoc Touch Imaging System and other digital imagers. A, western blot analysis for p44/42 MAPK (Erk1/2) expression was conducted using 2x serial dilutions (starting at 10 µg protein) of Jurkat cell lysate. The membranes were imaged on either the ChemiDoc Touch Imaging System or digital imagers from other vendors for a 15 sec exposure. As shown, the ChemiDoc Touch Imaging System is able to produce images with better definition and differentiation between closely spaced bands. B, the graph demonstrates the ability of the ChemiDoc Touch Imaging System to detect the same faint bands with greater intensity.

EASY, FLEXIBLE INTERACTION

Image Lab Touch Software takes the guesswork out of imaging and puts publication-quality images at your fingertips in seconds. Acquire images with a rapid 3-step workflow. Engage a full complement of digital tools to assess, select, and export your images.

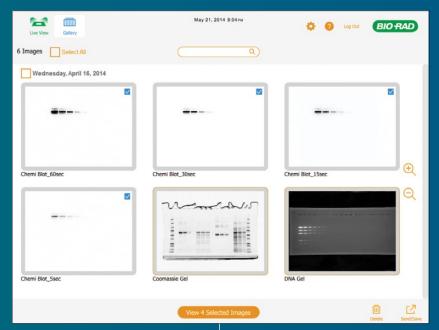




WESTERN BLOTTING CONSUMABLES

Assess and Export Images in the Gallery

The ChemiDoc Touch Imaging System has an intuitive interface to make reviewing, selecting, and exporting your images efficient and straightforward.



Gallery view enables you to peruse raw images



Pinch and zoom for a closer look



Compare up to 4 exposures side by side



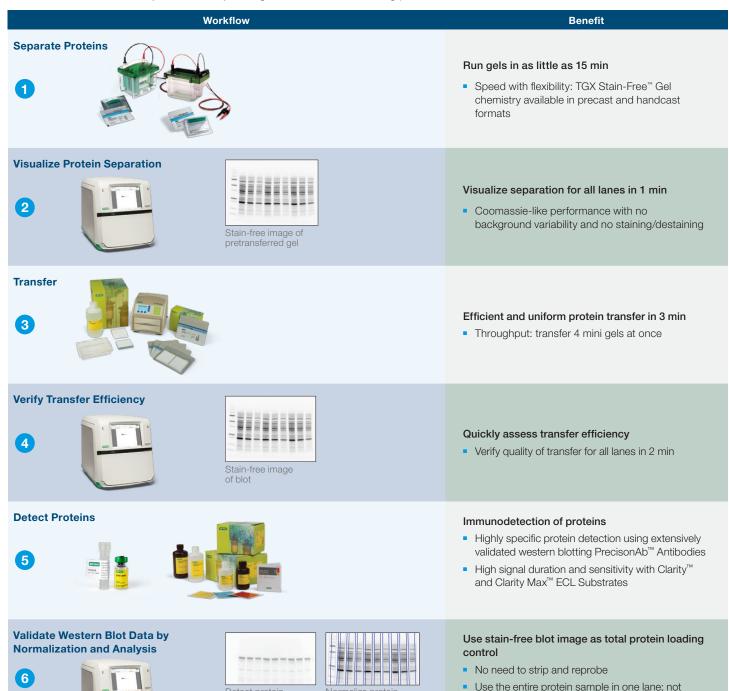
Export and print via USB or Ethernet connection

STAIN-FREE ENABLED

The ChemiDoc Touch Imaging System fully supports Bio-Rad's unique stain-free gel technology. Using the ChemiDoc Touch Imaging System as part of the V3 Western Workflow brings a new level of quality control and quantitation to the western blotting process, allowing multiple points at which to visualize, verify, and validate results.

V3 Western Workflow™

The V3 Western Workflow streamlines the western blotting protocol, incorporating stain-free in-gel chemistry to allow rapid fluorescent detection of proteins for gels and blots as well as the use of total protein normalization as a loading control. This improved workflow saves time and increases accuracy and reliability throughout the western blotting process.



Normalize protein

stain-free image of

blot from step 4

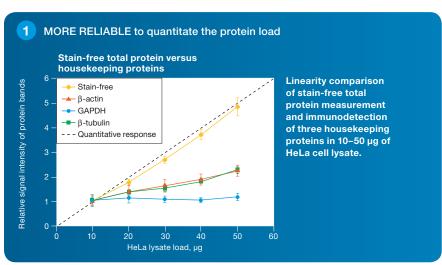
dependent on a single housekeeping protein

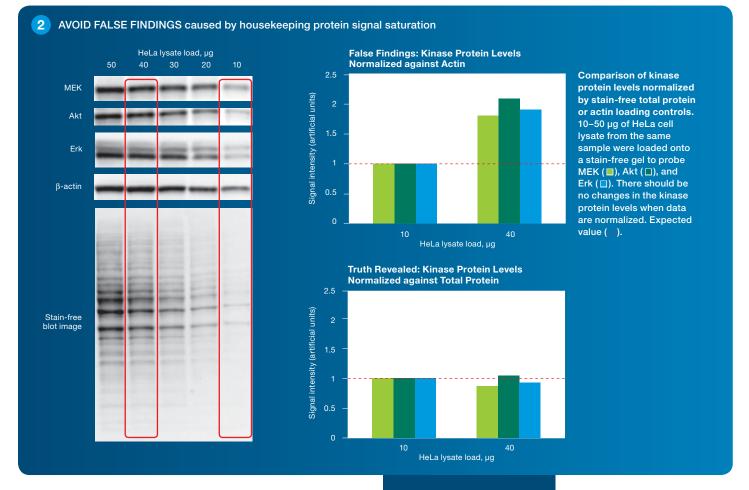
Reliable, accurate, and publishable quantitative data

Total Protein Normalization

Stain-free gel chemistry makes it possible to use total protein levels as a loading control rather than the housekeeping proteins used in traditional western blotting protocols. This negates the need to strip and reprobe the blot and avoids the attendant errors that can be introduced in this step.

Using total protein normalization produces a much greater linear dynamic range for measuring target protein levels. Housekeeping proteins such as β -actin, β -tubulin, or GAPDH are often very abundant in biological samples, which results in their signal being oversaturated compared to target proteins. Normalizing results to a total protein measurement corrects this problem, allowing a meaningful comparison even with low-abundance targets, and leads to far greater quantitative accuracy in measuring proteins of interest.

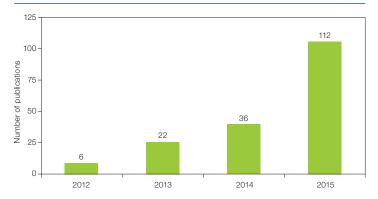




THE NEW GOLD STANDARD

Stain-free technology and, as a result, total protein normalization (TPN) are becoming increasingly popular in western blotting procedures. Major scientific journals are now asking researchers to conduct TPN preferentially over housekeeping normalization. In addition to the ChemiDoc Touch System, Bio-Rad also offers TGX Stain-Free Gels to support this more accurate quantitation method.

Peer-Reviewed Publications Using Total Protein Normalization



Stain-free total protein staining is a superior loading control to β -actin for western blots.

Gilda JE, Gomes AV (2013). Anal Biochem 440, 186-188.

V3 stain-free workflow for a practical, convenient, and reliable total protein loading control in Western blotting.

Posch A, Kohn J, Oh K, Hammond M, Liu N (2013). J Vis Exp, video ID 50948.

Rapid and precise engineering of the *Caenorhabditis elegans* genome with lethal mutation co-conversion and inactivation of NHEJ Repair.

Ward JD (2015). Genetics 199, 363-377.

Contractile properties and sarcoplasmic reticulum calcium content in type I and type II skeletal muscle fibres in active aged humans.

Lamboley CR, Wyckelsma VL, Dutka TL, McKenna MJ, Murphy RM, Lamb GD (2015). J Physiol 593, 2499–2514.

Alteration of mTOR signaling occurs early in the progression of Alzheimer disease (AD): Analysis of brain from subjects with pre-clinical AD, amnestic mild cognitive impairment and late-stage AD.

Tramutola A, Triplett JC, Di Domenico F, Niedowicz DM, Murphy MP, Coccia R, Perluigi M, Butterfield DA (2015). J Neurochem 133, 739–749.

Visit bio-rad.com/V3publications for a comprehensive list of publications using stain-free technology.

Incorporate TPN into Your Workflow with TGX Stain-Free Gel Chemistry

TGX Stain-Free products are designed with stain-free gel chemistry to provide superior gel performance and eliminate the need for staining. Compatible with standard sample and running buffers, TGX Stain-Free products allow you to

- Quickly visualize proteins no staining required
- Run gels in as little as 15 min
- Efficiently transfer proteins in as little as 3 min

They are available in multiple formats for your convenience:

 TGX Stain-Free Precast Gels — Mini-PROTEAN[®] (mini) and Criterion[™] (midi) Gels optimized for western blotting



Mini-PROTEAN Gels — bio-rad.com/MiniStainFree1
Criterion Gels — bio-rad.com/MidiStainFree1

 TGX Stain-Free[™] FastCast[™] Acrylamide Solutions — premixed acrylamide solutions for hand casting polyacrylamide gels



 ${\sf FastCast\ Solutions-bio-rad.com/SFFastCast1}$

WESTERN BLOTTING CONSUMABLES

The ChemiDoc Touch Imaging System is part of Bio-Rad's range of products to improve the entire western blot process, from immunoprecipitation all the way through to data analysis. These consumables provide workflow optimizations and better results for a variety of laboratory needs.

Immunoprecipitation with SureBeads[™] Magnetic Beads

For protein complex pull-down and isolation of low-abundance targets

- Faster and easier way to immunoprecipitate say yes to magnetization, no to centrifugation
- Patented surface chemistry enables proper antibody orientation, which maximizes antigen binding capacity
- Ergonomically designed 16-tube SureBeads Magnetic Rack has strong separable magnets to minimize sample handling and is fast, easy to use, and affordable.

bio-rad.com/MagneticIP



Precision Plus Protein[™] Standards

Designed for accurate molecular weight estimation

- Precision Plus Protein Dual Color Standards brighter for easier target protein identification and can yield stronger band intensity after blot processing
- Precision Plus Protein All Blue Standards recommended for colorimetric visualization on stain-free gels
- Precision Plus Protein Unstained Standards provide stain-free visible ladder

bio-rad.com/PrecisionPlus1



Prepacked Transfer Consumables

All the resources needed for a fast and efficient transfer process

- Trans-Blot[®] Turbo[™] Transfer Packs reduce setup time to 1 min from the opening of the gel cassette to the start of your transfer
- Ready-to-assemble (RTA) transfer kits provide all consumables needed to transfer 40 blots, including transfer buffer and transfer stacks available with nitrocellulose, PVDF, or low fluorescence PVDF membranes. A cost-effective and easy-to-use solution for protein transfers

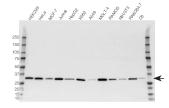
bio-rad.com/TransBlotTurbo

PrecisionAb Antibodies

Rigorously validated to work every time

- Highly specific and extensively validated with up to 12 cell lysates
- Strict quality control for lot-to-lot consistency
- Trial size for cost effective evaluation
- Positive control lysate to facilitate target identification

bio-rad.com/WBantibodies



Clarity and Clarity Max Western ECL Substrates

Compatible with any HRP-conjugated secondary antibody

The perfect solution for detecting high- and low-expressing proteins, even when taking multiple exposures.

- Clarity Substrate moderate sensitivity, long signal duration, and two-year shelf life at room temperature
- Clarity Max Substrate for when you need high sensitivity

bio-rad.com/ClarityECL



Specifications

Specifications	
Automation Capabilities	
Smart Tray Technology	ChemiDoc Touch Imaging System automatically recognizes your application-specific tray and adjusts imaging parameters and software options accordingly
Autofocus	Precalibrated focus for any zoom setting or sample height
Auto-exposure	two user-defined modes (rapid or optimal auto-exposure) for chemiluminescence two user-defined modes (faint or intense bands) for
	nonchemiluminescence applications
Image flat fielding	Dynamic; precalibrated and optimized for every application
Hardware Specifications	;
Touch screen functionality	Multitouch capable (4 points) 12.1" display
Maximum image area (L x W)	16.8 x 21 cm
Illumination source	Trans-UV, 302 nm (standard)
	Epi-white (standard)
	Trans-white (optional)
	Trans-blue (optional)
Detector	Cooled CCD, 6 megapixels
Camera cooling temperature	-25°C
Filter holder	Two positions (one for standard filter, one without filter for chemiluminescence)
Emission filter	Standard filter to perform protein and DNA gel and blot imaging
Dynamic range	>4 orders of magnitude
Data output	16-bit or 8-bit; SCN, TIFF, JPEG image files
Instrument weight	35 kg (78 lbs)
Instrument size (L x W x H)	61 x 51 x 53 cm
Operating voltage	100-250 V
Operating temperature	10-28°C

10-85% relative humidity (noncondensing)

Ordering Information

1707581 Software 1709690

Catalog #	Description
1708370	ChemiDoc Touch Imaging System, includes internal computer, 12" touch-screen display, camera, Image Lab Touch Software, chemi/UV/stain-free sample tray, Clarity Western ECL Substrate, Precision Plus Protein Dual Color Standards
1708381	ChemiDoc Touch V3 Western Workflow for Mini Gels, includes ChemiDoc Touch Imaging System with Image Lab Touch Software, chemi/UV/stain-free sample tray, 50 Any kD™ Mini-PROTEAN TGX Stain-Free Precast Gels, SDS-PAGE accessories, Clarity Western ECL Substrate, Precision Plus Protein Dual Color Standards, Mini-PROTEAN Tetra Cell, Trans-Blot® Turbo™ Transfer Starter System, 50 PVDF transfer packs for mini gels
1708382	ChemiDoc Touch V3 Western Workflow for Midi Gels, includes ChemiDoc Touch Imaging System with Image Lab Touch Software, chemi/UV/stain-free sample tray, 50 4–20% Criterion TGX Stain-Free Precast Gels, SDS-PAGE accessories, Clarity Western ECL Substrate, Precision Plus Protein Dual Color Standards, Criterion Cell, Trans-Blot Turbo Transfer Starter System, 50 PVDF transfer packs for midi gels
Accessories	
1708372	White sample tray, for gels stained with Coomassie Blue, copper, silver, or zinc stains
1708373	Blue sample tray, with viewing goggles, for gels stained with GelGreen or any SYBR® Stains
1708373 1708374	
	GelGreen or any SYBR® Stains Chemi/UV/stain-free sample tray, for chemiluminescent blots, stain-free gels/blots, and gels stained with ethidium bromide,
1708374	GelGreen or any SYBR® Stains Chemi/UV/stain-free sample tray, for chemiluminescent blots, stain-free gels/blots, and gels stained with ethidium bromide, SYPRO Ruby, Oriole™, GelRed, and SYBR® Stains. UV safety shield, to protect against UV light exposure during
1708374 1708375	GelGreen or any SYBR® Stains Chemi/UV/stain-free sample tray, for chemiluminescent blots, stain-free gels/blots, and gels stained with ethidium bromide, SYPRO Ruby, Oriole™, GelRed, and SYBR® Stains. UV safety shield, to protect against UV light exposure during band excision Gel alignment templates, for consistent placement of gels
1708374 1708375 1708376	GelGreen or any SYBR® Stains Chemi/UV/stain-free sample tray, for chemiluminescent blots, stain-free gels/blots, and gels stained with ethidium bromide, SYPRO Ruby, Oriole™, GelRed, and SYBR® Stains. UV safety shield, to protect against UV light exposure during band excision Gel alignment templates, for consistent placement of gels and blots Holder for sample trays and UV shield ChemiDoc Touch IQ/OQ protocols, for installation qualification/
1708374 1708375 1708376 1708377 1708378	GelGreen or any SYBR® Stains Chemi/UV/stain-free sample tray, for chemiluminescent blots, stain-free gels/blots, and gels stained with ethidium bromide, SYPRO Ruby, Oriole™, GelRed, and SYBR® Stains. UV safety shield, to protect against UV light exposure during band excision Gel alignment templates, for consistent placement of gels and blots Holder for sample trays and UV shield ChemiDoc Touch IQ/OQ protocols, for installation qualification/operational qualification
1708374 1708375 1708376 1708377 1708378 1708097	GelGreen or any SYBR® Stains Chemi/UV/stain-free sample tray, for chemiluminescent blots, stain-free gels/blots, and gels stained with ethidium bromide, SYPRO Ruby, Oriole™, GelRed, and SYBR® Stains. UV safety shield, to protect against UV light exposure during band excision Gel alignment templates, for consistent placement of gels and blots Holder for sample trays and UV shield ChemiDoc Touch IQ/OQ protocols, for installation qualification/ operational qualification Standard 302 nm UV lamps, pkg of 6
1708374 1708375 1708376 1708377 1708378	GelGreen or any SYBR® Stains Chemi/UV/stain-free sample tray, for chemiluminescent blots, stain-free gels/blots, and gels stained with ethidium bromide, SYPRO Ruby, Oriole™, GelRed, and SYBR® Stains. UV safety shield, to protect against UV light exposure during band excision Gel alignment templates, for consistent placement of gels and blots Holder for sample trays and UV shield ChemiDoc Touch IQ/OQ protocols, for installation qualification/operational qualification

Visit bio-rad.com/CDTinfo to learn more about the ChemiDoc Touch Imaging System.

for viewing images and 1-D analysis

SYPRO and SYBR are trademarks of Life Technologies Corporation. GelGreen and GelRed are trademarks of Biotium, Inc. Mitsubishi is a trademark of Mitsubishi Companies. Mac is a trademark of Apple Inc.

Image Lab Software, stand-alone version, PC or Mac,

Bio-Rad Laboratories, Inc. is licensed by Life Technologies Corporation to sell SYPRO products for research use only under U.S. Patent Number 5,616,502.

Precision Plus Protein Standards are sold under license from Life Technologies Corporation, Carlsbad, CA for use only by the buyer of the product. The buyer is not authorized to sell or resell this product or its components.



Operating humidity

Bio-Rad Laboratories, Inc.

Life Science Group

Web site bio-rad.com USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 43 1 877 89 01 177 Belgium 32 (0)3 710 53 00 Brazil 55 11 3065 7550 Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 45 44 52 10 00 Finland 358 09 804 22 00 France 33 01 47 95 69 65 Germany 49 89 31 884 0 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300

Israel 972 03 963 6050 Italy 39 02 216091 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 31 (0)318 540 666

New Zealand 64 9 415 2280 Norway 47 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04

Singapore 65 6415 3188 South Africa 27 (0) 861 246 723 Spain 34 91 590 5200 Sweden 46 08 555 12700 Switzerland 41 026 674 55 05

Taiwan 886 2 2578 7189 Thailand 66 662 651 8311 United Arab Emirates 971 4 8187300 United Kingdom 44 020 8328 2000

Bulletin 6517 Ver D US/EG 16-0522 0516 Sig 1215

