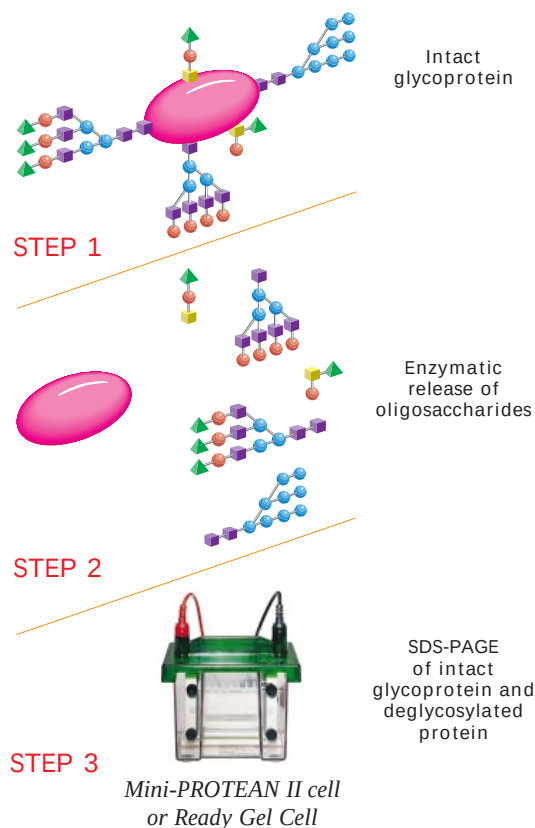


KIT 2⁺

HOW DO I REMOVE OLIGOSACCHARIDES FROM MY GLYCOPROTEIN?

The Enzymatic Deglycosylation and Deglycosylation Enhancement kits provide a gentle method that leaves both carbohydrate and peptide intact for further protein and carbohydrate analysis. The Enzymatic Deglycosylation Kit contains PNGase F, NANase II, and O-Glycosidase, which remove all N-linked oligosaccharides and the sialylated and non-sialylated O-linked core structure of Gal(β-1,3)GalNAc(α1). The Enhancement Kit, which contains GALase III and HEXase I, is used in combination with the Deglycosylation Kit to remove additional O-linked core structures containing β1,4 linked galactose and N-acetylglucosamine.



SPECIFICATIONS

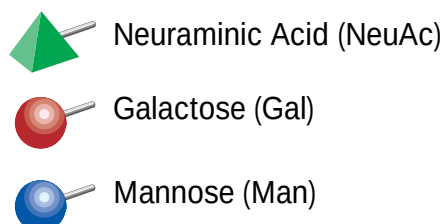
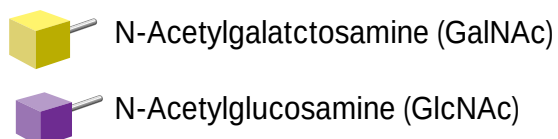
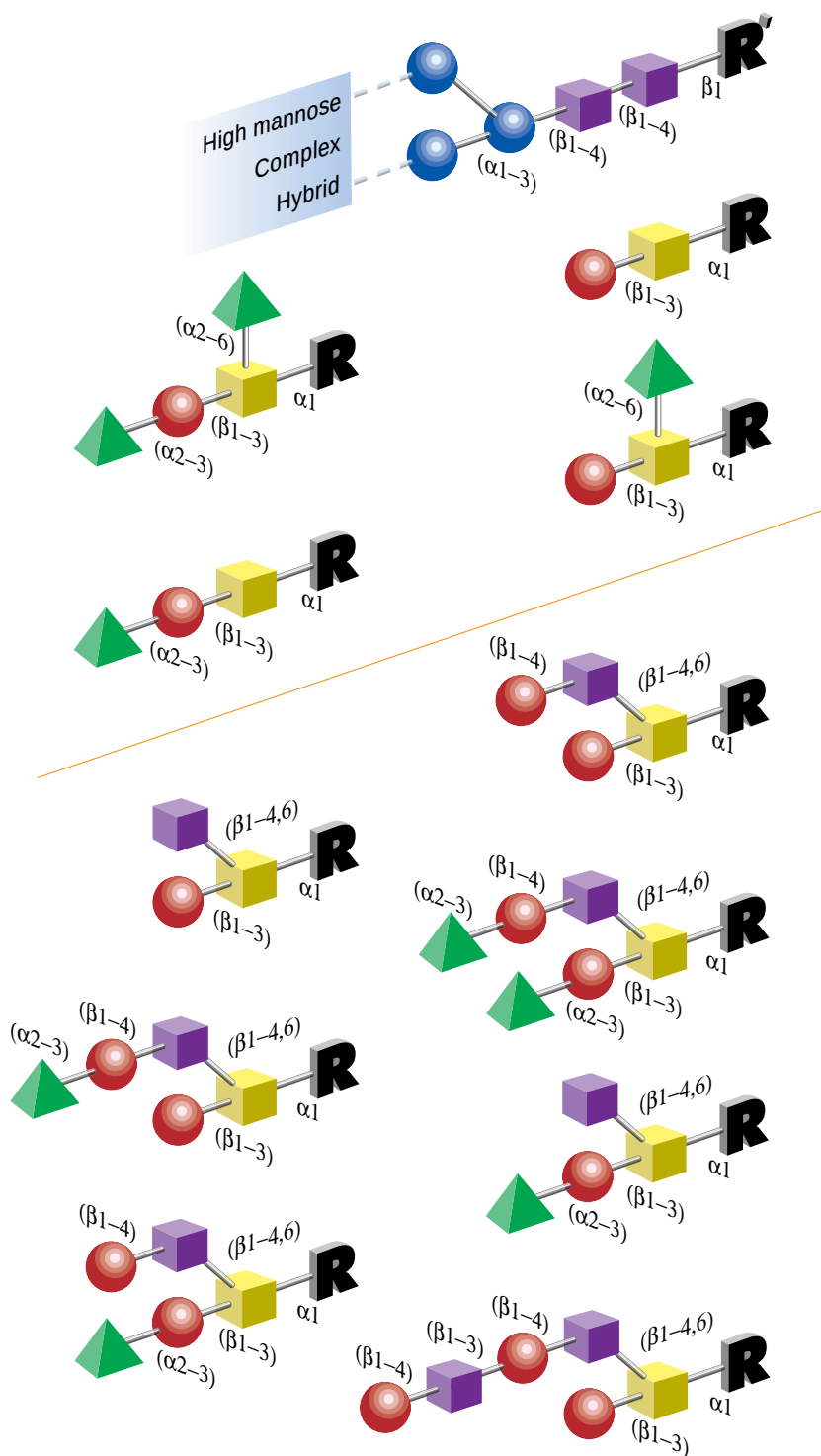
ENZYME	CONCENTRATION	SPECIFICITY
Enzymatic Deglycosylation Kit		
NANase II	5 U/ml	All α(2-3, 6) N-acetylneuraminic acids
O-Glycosidase DS	1 U/ml	All Ser/Thr linked Gal (β1-3)GalNAc(α1)
PNGase F	2.5 U/ml	All Asn-linked oligosaccharides
Deglycosylation Enhancement Kit		
GALase III	1.5 U/ml	Non-reducing terminal (β1-4) linked galactose from complex carbohydrates
HEXase I	8 U/ml	Non-reducing terminal (β1-2, 3, 4, 6) N-acetyl-glucosamine from complex carbohydrates

ORDERING INFORMATION

CATALOG NUMBER	PRODUCT DESCRIPTION
170-6500	Enzymatic Deglycosylation Kit
170-6508	Deglycosylation Enhancement Kit
170-6880	HEXase I, 42 U/ml, 40 µl
170-6881	O-Glycosidase DS, 1 U/ml, 40 µl
170-6882	NANase II, 5 U/ml, 40 µl
170-6883	PNGase F, 2.5 U/ml, 40 µl
170-6513	GALase III, 1.5 U/ml, 40 µl

N- and O-Linked Cores Removed with the Enzymatic Deglycosylation Kit

Additional O-Linked Cores Removed with the Deglycosylation Enhancement Kit in Combination with the Enzymatic Deglycosylation Kit



BIO-RAD

**Bio-Rad
Laboratories**

Life Science
Group

U.S. (800) 4BIORAD • California (510) 741-1000 • Australia 02-9914-2800 • Austria (1)-877 89 01 • Belgium 09-385 55 11 • Canada (905) 712-2771 • China (01) 2046622 • Denmark 39 17 9947 • Finland 90 804 2200 • France (1) 49 60 68 34 • Germany 089 318 84-0 • India 91-11-461-0103 • Italy 02-21609 1 • Japan 03-5811-6270 • Hong Kong 7893300 • The Netherlands 0318-540666 • New Zealand 09-443 3099 • Singapore (65) 272-9877 • Spain (91) 661 70 85 • Sweden 46 (0) 8-735 83 00 • Switzerland 01-809 55 55 • United Kingdom 0800 181134