



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

SG-Chymotrypsin™

(Cat. # 786-13)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

SG-Chymotrypsin™ is a sequencing grade serine endopeptidase, which predominantly cleaves peptide bonds on the carboxy side of tyrosine, phenylalanine and tryptophan. In addition, it has a low catalytic activity against the carboxy side of leucine, methionine, alanine, aspartic and glutamic acids, although at a much lower rate. It is therefore recommended to always use the shortest digestion time possible. SG-Chymotrypsin™ is first treated with TLCK and then subjected to an extensive purification process to remove contaminating protease and chymotryptic autolysis by-products. The highly purified enzyme is then chemically modified to increase its resistance to autolysis and increase its stability. The modified enzyme retains >80% of its activity after 6 hours incubation at 30°C in reaction buffer and >70% of activity after 24 hours incubation under the same conditions. The chemically modified SG-Chymotrypsin™ is stable in denaturing agents (see Table) and therefore can be used to digest difficult to solubilize proteins.

Denaturing Agent	Concentration	% Enzyme Activity Retained
Control	-	100
Urea	0.50M	100
	1.00M	100
	2.00M	100
	3.00M	100
	4.00M	100
Guanidine-HCl	0.05M	100
	0.10M	100
	0.25M	100
	0.50M	11

ITEM(S) SUPPLIED (CAT. # 786-13)

Description	Size
SG-Chymotrypsin™	2 vials, 5µg/vial
Digestion Buffer (CHY)	2ml

STORAGE CONDITIONS

It is shipped at ambient temperature. Upon arrival, store at -20°C and is stable for 1 year.

PREPARATION BEFORE USE

NOTE: SG-Chymotrypsin TM is supplied lyophilized, 5 µg/vial.

Reconstitute the enzyme with 100 µl sterile water to produce a concentration of 100 ng/µl. Combine 5 µl of 100 ng/µl solution with 95 µl Digestion Buffer (CHY) to yield 5 ng/µl chymotrypsin. Reconstituted enzyme is stable for 1 month at -20 °C, repeated freeze thawing is not recommended.

NOTE: If reduction and alkylation are needed 250 mM dithiothreitol and 625 mM iodoacetamide are recommended.

PROTOCOL

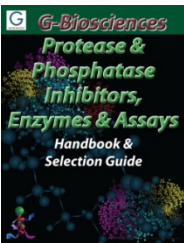
For optimal digestion make sure protein sample is either prepared or equilibrated in 50 mM Tris-HCl, pH 8.0, 0.1 mM CaCl₂.

1. For protein fragmentation, 5 ng/µl SG-Chymotrypsin TM is typically added to the protein at a ratio of 1:20 to 1:100 enzyme to protein, by weight.
2. The incubation is allowed to proceed at 37 °C for 2-10 hours, but can be extended to 24 hours in some applications.

NOTE: It is recommended to choose a ratio of enzyme to protein to allow for the shortest incubation time possible, to reduce or eliminate the catalyzed hydrolysis of peptide bonds with non-aromatic amino acid residues.

RELATED PRODUCTS

Download our Protease & Phosphatase Inhibitors, Enzyme & Assays Handbook.



<http://info.gbiosciences.com/protease-phosphatase-inhibitors-enzymes-assay-handbook>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 5/6/2016 1A



www.GBiosciences.com