

Certificate of Analysis

IdeZ Protease, Frozen, 2,000 units

Cat.# V8342

Includes:

Part#	Name	Size
V834B	IdeZ Protease, Frozen	2,000 units

Description: IdeZ (Immunoglobulin-degrading enzyme from *Streptococcus equi* subspecies *zooepidemicus*) is an engineered recombinant protease overexpressed in *Escherichia coli*. IdeZ Protease specifically cleaves IgG molecules below the hinge region to yield F(ab)₂ and Fc fragments.

Biological Source: Recombinant strain of *E. coli*.

Molecular Weight: IdeZ Protease has a molecular weight of approximately 38kDa.

Storage Conditions: Store product below -10°C. After initial thawing, IdeZ Protease may be stored at 2-10°C for up to 60 days.

Concentration: 50 units/μl in 50mM sodium phosphate, 150mM NaCl (pH 6.6).

Expiration Date: See product label for expiration date.

Unit Definition: One unit will cleave ≥95% of 1μg of recombinant monoclonal IgG in 30 minutes at 37°C.

Part# 9PIV8342

Revised 2/17



AF9PIV83420217V8342



Promega

Promega Corporation

2800 Woods Hollow Road	
Madison, WI 53711-5399	USA
Telephone	608-274-4330
Toll Free	800-356-9526
Fax	608-277-2516
Internet	www.promega.com

Quality Control Assays

This lot passes the following Quality Control specifications:

Purity: ≥95% as determined by SDS-PAGE analysis.

Activity: Unit activity for each lot of IdeZ Protease is confirmed by the cleavage of recombinant monoclonal IgG as analyzed by SDS-PAGE.

PRODUCT USE LIMITATIONS, WARRANTY DISCLAIMER

Promega manufactures products for a number of intended uses. Please refer to the product label for the intended use statements for specific products. Promega products contain chemicals which may be harmful if mis-used. Due care should be exercised with all Promega products to prevent direct human contact.

Each Promega product is shipped with documentation stating specifications and other technical information. Promega products are warranted to meet or exceed the stated specifications. Promega's sole obligation and the customer's sole remedy is limited to replacement of products free of charge in the event products fail to perform as warranted. Promega makes no other warranty of any kind whatsoever, and SPECIFICALLY DISCLAIMS AND EXCLUDES ALL OTHER WARRANTIES OF ANY KIND OR NATURE WHATSOEVER, DIRECTLY OR INDIRECTLY, EXPRESS OR IMPLIED, INCLUDING, WITHOUT LIMITATION, AS TO THE SUITABILITY, PRODUCTIVITY, DURABILITY, FITNESS FOR A PARTICULAR PURPOSE OR USE, MERCHANTABILITY, CONDITION, OR ANY OTHER MATTER WITH RESPECT TO PROMEGA PRODUCTS. In no event shall Promega be liable for claims for any other damages, whether direct, incidental, foreseeable, consequential, or special (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including negligence) or strict liability arising in connection with the sale or the failure of Promega products to perform in accordance with the stated specifications.

Signed by:

R. Wheeler, Quality Assurance

© 2015-2017 Promega Corporation. All Rights Reserved.

ISOQUANT is a registered trademark of Promega Corporation. Magne is a trademark of Promega Corporation.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All prices and specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.

Protocol for Antibody Fragmentation

1. Add the desired amount of IgG (up to 5mg) in digestion buffer or other compatible buffer¹.
2. Add IdeZ Protease to the IgG sample:
 - Add 1 unit of IdeZ Protease per 1µg of IgG to be digested.
 - For example, add 1µl (50 units) of IdeZ Protease to digest 50µg of IgG.
3. Incubate sample at 37°C for 30–60 minutes².

¹IdeZ Protease is most active in buffers at or near neutral pH. The recommended digestion buffer is 50mM sodium phosphate, 150mM NaCl (pH 6.6), but many common biological buffers such as Tris or PBS also can be used. Buffers outside this pH range (e.g., acetate buffers) may be used, but the incubation time and/or enzyme amount may have to be optimized on a case-by-case basis.

²Mouse IgG2a and IgG3 typically require 2–4 hours for complete digestion. In some cases, human IgG2 may require incubation beyond 60 minutes.

Notes

- IgG concentration ideally should be in the range of 1–20mg/ml.
- IdeZ Protease efficiently cleaves human, humanized, chimeric, sheep, rabbit and monkey IgGs as well as mouse IgG2a and IgG3. IdeZ Protease also will cleave many Fc-fusion proteins as well as antibody-drug conjugates (ADCs).
- IdeZ Protease does not cleave mouse IgG1/IgG2b, rat, porcine, bovine or goat IgG. IdeZ Protease also does not cleave non-IgG isotypes including IgA, IgM, IgD and IgE.
- IdeZ Protease cleaves at a single site below the hinge at the same location as IdeS Protease (Cat.# V7511, V7515).
- The IdeZ Protease has a histidine tag for easy removal.
- For downstream analysis of fragments by LC-MS, best results are obtained following reduction and denaturing of the IdeZ-digested fragments.
- Purified F(ab)₂ fragments can be obtained after digestion by incubating the digest for 30–60 minutes with Magne™ Protein A Beads (Cat.# G8781).
- IdeZ Protease can be used in the same reaction as PNGase F (Cat.# V4831) to perform fragmentation and removal of Fc glycans in a single step using the recommended digest buffer. Ten units of PNGase F may be sufficient to remove Fc glycans from 50µg of IgG in 2 hours at 37°C, although some optimization of PNGase F amount and incubation time may be required.

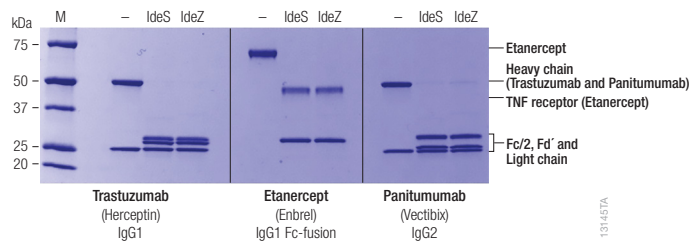


Figure 1. Digestion of panel of therapeutic IgGs and Fc-fusion proteins. Therapeutic IgG or fusion protein (50µg) was digested with 50 units of IdeS Protease or IdeZ Protease for 30 minutes at 37°C in a final volume of 25µl in the recommended digestion buffer (50mM sodium phosphate, 150mM NaCl [pH 6.6]). Undigested controls and samples were analyzed by SDS-PAGE under reducing conditions.

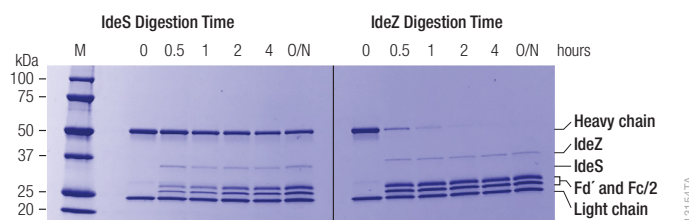


Figure 2. Digestion of recombinant mouse IgG2a with IdeS Protease and IdeZ Protease. Recombinant mIgG2a anti-hCD20 (50µg; Invivogen) was digested with 50 units of IdeS Protease or IdeZ Protease for various time points at 37°C in a final volume of 25µl in the recommended digestion buffer (50mM sodium phosphate, 150mM NaCl [pH 6.6]). IgG2a is completely digested in 2 hours with IdeZ Protease but is only partially digested with IdeS Protease even after an overnight (O/N) digestion.

Related Products

Product	Size	Cat.#
IdeZ Protease (lyophilized)	5,000 units	V8341
	25,000 units	
	(5 × 5,000 units)	V8345
IdeS Protease (lyophilized)	5,000 units	V7511
	25,000 units	
	(5 × 5,000 units)	V7515
IdeS Protease, Frozen	2,000 units	V7512
PNGase F	500 units	V4831
Magne™ Protein A Beads, 20% Slurry	1ml	G8781
ISOQUANT® Isoaspartate Detection Kit*	100 assays	MA1010
Protein Deglycosylation Mix	20 reactions	V4931
Trypsin/Lys-C Mix, Mass Spec Grade	100µg	V5072
Trypsin Gold, Mass Spectrometry Grade	100µg	V5280
Sequencing Grade Modified Trypsin	100µg (5 × 20µg)	V5111
Sequencing Grade Modified Trypsin, Frozen	100µg (5 × 20µg)	V5113
rLys-C, Mass Spec Grade	15µg	V1671
Asp-N, Sequencing Grade	2µg	V1621
Glu-C, Sequencing Grade	50µg (5 × 10µg)	V1651

*Not For Medical Diagnostic Use.