

Data Sheet

IDE

Human, recombinant, N-terminal His-tag

Catalog #: 70002

Lot #: 140805

Conc.: 0.84 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 3 mM DTT, and 105 mM imidazole

Stability: >6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted protein is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. Farris, W., et al., *Proc. Natl. Acad. Sci. (USA)*. 2003; **100(7)**: 4162-4167.
2. Qiu, W.Q., et al., *J. Biol. Chem.* 1998; **273**: 32730-32738.

Description: Human insulin-degrading enzyme (IDE), also known as insulysin, Abeta-degrading protease, insulin protease, and insulinase, GenBank Accession No. NM_004969, a.a. 42-1019(end) with N-terminal His-tag, expressed in a Baculovirus Sf9 cell expression system. MW = 114 kDa.

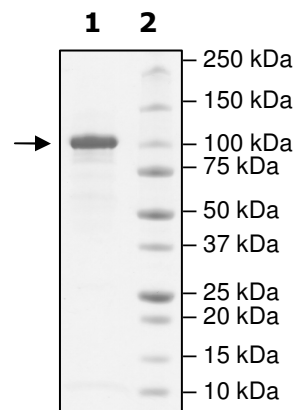
Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
3.7 μg IDE
Lane 2:
Protein Marker

Purity: $\geq 80\%$
MW: 114 kDa



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com