

## Data Sheet

### HEXB

Human, recombinant, C-terminal FLAG-His-tags

**Catalog #:** 71088

Lot#: 130503-H

Conc.: 0.23 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 150 mM imidazole and 20% glycerol**Stability:** >6 months at -80 °C**References:**

1. Mahuran, D.J., *et al.*, *Biochim Biophys Acta*. 1991 Feb 22;**1096(2)**:87-94.
2. Mark, B.L., *et al.*, *J Mol Biol*. 2003;**327**:1093-1109.

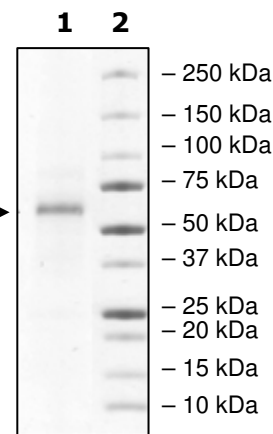
**Description:**

Human hexosaminidase B (beta polypeptide), also known as HEXB, (GenBank Accession No. NM\_000521), full length (a.a. 1-556(end)) with C-terminal FLAG-His-tags, MW = 64.9 kDa, expressed in a baculovirus infected Sf9 cell expression system.

**Application:**

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

## Quality Assurance

**4-20% SDS-PAGE  
Coomassie staining****Lane 1:**  
2.8 µg HEXB →  
**Lane 2:**  
Protein Marker**MW:** 64.9 kDa  
**Purity:** ≥82%

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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court West, Suite E  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Sequence:**

MELCGLGLPRPPMLLALLLATLLAAMLALLTQVALVVQVAEAAARAPSVSAKPGPALWP  
LPLSVKMPNLLHLAPENFYISHSPNSTAGPSCTLLEAFRRYHGYIFGFYKWHHEPAE  
FQAKTQVQQLLVSITLQSECDAFPNISSDESYLLVKEPVAVLKANRVWGALRGLETFS  
QLVYQDSYGTFTINESTIIDSPRFSHRGILIDTSRHYLPVKIILKTLDAMAFNKFNVLHWHIV  
DDQSFPYQSITFPELSNKGSYLSHVYTPNDVRMVIEYARLRGIRVLPEFDTPGHTLSWG  
KGQKDLLTPCYSRQNKLDSEFGPINPTLNTTYSFLTTFKEISEVFPDQFIHLGGDEVEFK  
CWESNPKIQDFMRQKGFGTDFKKLESFYIQKVLDIATINKGSIVWQEVFDDKAKLAPGTI  
VEVWKDSAYPEELSRVTASGFPVILSAPWYLDLISYGQDWRKYYKVEPLDFGGTQKQK  
QLFIGGEACLWGEYVDATNLTPRLWPRASAVGERLWSSKDVRDMDDAYDRLTRHRCR  
MVERGIAAQPLYAGYCNHENM**DYKDDDDKHHHHH**

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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)