

PRODUCT INFORMATION



GBA3 (human, recombinant)

Item No. 33991

Overview and Properties

Synonyms: cBGL1, Cytosolic β -Glucosidase, Cytosolic β -Glucosidase-like Protein-1, Klotho-related Protein, KLRP

Source: Active recombinant human C-terminal His-tagged GBA3 expressed in insect cells

Amino Acids: 1-469 (full length)

Uniprot No.: Q9H227

Molecular Weight: 55 kDa

Storage: -80°C (as supplied)

Stability: ≥ 1 year

Purity: $\geq 90\%$ estimated by SDS-PAGE

Supplied in: Lyophilized from sterile 20 mM Tris, 300 mM sodium chloride, pH 7.5, with 10% glycerol, 5% trehalose, 5% mannitol, and 0.01% tween-80

Endotoxin Testing: < 1.0 EU/ μ g, determined by the LAL endotoxin assay

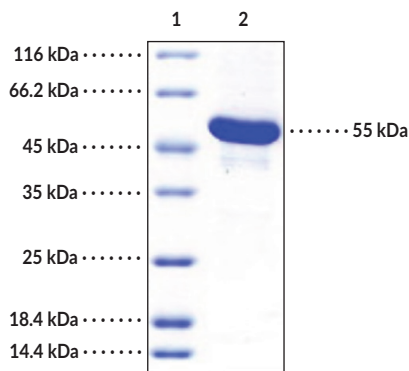
Protein Concentration: *batch specific* mg/ml

Activity: Measured by its ability to hydrolyze 4-methylumbelliferyl- β -D glucopyranoside.

Specific Activity: *batch specific* U/mg

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



Lane 1: MW Markers
Lane 2: GBA3

SDS-PAGE Analysis of GBA3. This protein has a calculated molecular weight of 55 kDa.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/14/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Cytosolic β -glucosidase (GBA3) is a neutral β -glucocerebrosidase.^{1,2} It is composed of a single glycoside hydrolase domain and hydrolyzes glucosyl- and galactosylceramides. GBA3 is expressed in the liver, small intestine, colon, spleen, and kidneys.³ *In vitro*, knockdown of GBA3 inhibits mitotic arrest and promotes resistance to paclitaxel (Item No. 10461) in colon cancer cells.⁴ Tumor levels of GBA3 are lower in patients with hepatocellular carcinoma and decreased expression is associated with poor prognosis.⁵ Unlike GBA1, polymorphisms in GBA3 are not associated with Gaucher disease, a lysosomal storage disorder.⁶ Cayman's GBA3 (human, recombinant) protein can be used for enzyme activity assays. This protein consists of 480 amino acids and has a calculated molecular weight of 55 kDa.

References

1. Hayashi, Y. and Ito, M. Klotho-related protein KLRP: Structure and functions. *Vitamins and Hormones* **101**, 1-16 (2016).
2. Astudillo, L., Therville, N., Colacios, C., *et al.* Glucosylceramidases and malignancies in mammals. *Biochimie* **125**, 267-280 (2016).
3. Yahata, K., Mori, K., Arai, H., *et al.* Molecular cloning and expression of a novel klotho-related protein. *J. Mol. Med. (Berl.)* **78(7)**, 389-394 (2000).
4. Swanton, C., Marani, M., Pardo, O., *et al.* Regulators of mitotic arrest and ceramide metabolism are determinants of sensitivity to paclitaxel and other chemotherapeutic drugs. *Cancer Cell* **11(6)**, 498-512 (2007).
5. Ying, J.F., Zhang, Y.N., Song, S.S., *et al.* Decreased expression of GBA3 correlates with a poor prognosis in hepatocellular carcinoma patients. *Neoplasma* **67(5)**, 1139-1145 (2020).
6. Beutler, E., Beutler, L., and West, C. Mutations in the gene encoding cytosolic β -glucosidase in Gaucher disease. *J. Lab. Clin. Med.* **144(2)**, 65-68 (2004).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM