

PRODUCT INFORMATION



BTBD11 (human, recombinant)

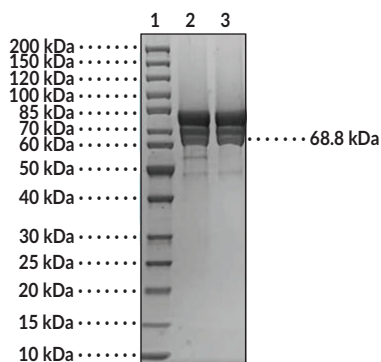
Item No. 45175

Overview and Properties

Synonyms:	ABTB3, Ankyrin Repeat- and BTB/POZ Domain-containing Protein 3, BTB/POZ Domain-containing Protein 11
Source:	Recombinant human N-terminal His-Strep II-tagged BTBD11 expressed in insect cells
Amino Acids:	2-1,104 (full length)
Uniprot No.:	A6QL63
Molecular Weight:	68.8 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	≥90% estimated by SDS-PAGE
Supplied in:	50 mM Tris-HCl, pH 7.5, with 200 mM sodium chloride, 20% glycerol, and 1 mM DTT
Endotoxin Testing:	<1.0 EU/μg, determined by the LAL endotoxin assay
Protein Concentration:	<i>batch specific</i> mg/ml

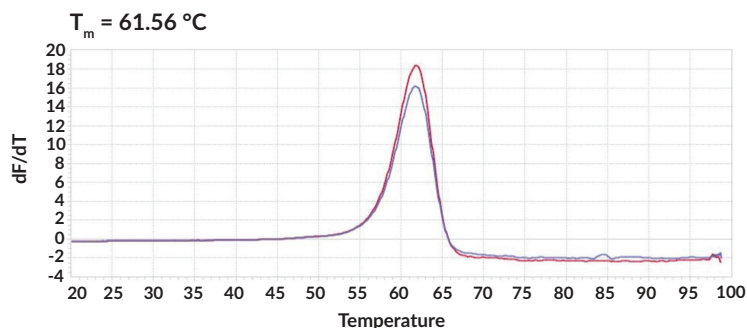
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: BTBD11 (2 μg) (reduced)
Lane 3: BTBD11 (2 μg) (non-reduced)

SDS-PAGE Analysis of BTBD11. This protein has a calculated molecular weight of 68.8 kDa.



Protein Melting Thermal Profile of BTBD11.

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.

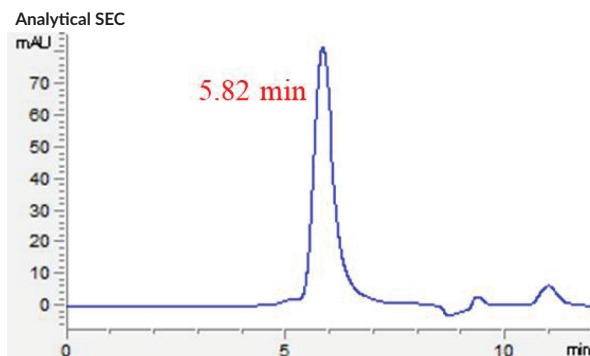
WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/18/2026

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



Analytical size exclusion chromatography of protein BTBD11.

Description

BTB/POZ domain-containing protein 11 (BTBD11) is a BTB-domain containing protein.¹ It is composed of a disordered region, five ankyrin repeats, a BTB domain, and a PDZ-binding motif (PBM). Btbd11 is expressed in cortical and hippocampal inhibitory interneurons. It binds to postsynaptic density protein 95 (Psd-95) and is involved in regulation of glutamatergic synapses and neuronal circuit function. Knockout of the gene encoding Btbd11, *Abtb3*, in inhibitory neurons expressing the vesicular inhibitory amino acid transporter (VIAAT), also known as VGAT, increases rearing events in mice and decreases anxiety-like behavior in male mice. A SNP in *ABTB3* is associated with risk of human papillomavirus-negative (HPV-negative) oropharyngeal cancer.² Cayman's BTBD11 (human, recombinant) protein has a calculated molecular weight of 68.8 kDa.

References

1. Bygrave, A.M., Sengupta, A., Jackert, E.P., *et al.* Btbd11 supports cell-type-specific synaptic function. *Cell Rep.* **42**(6), 112591 (2023).
2. Ferreiro-Iglesias, A., McKay, J.D., Brenner, N., *et al.* Germline determinants of humoral immune response to HPV-16 protect against oropharyngeal cancer. *Nat. Commun.* **12**(1), 5945 (2021).

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