

HSD17B10 Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog # H23-30G

Lot # U316-2

Product Description

Recombinant full-length human HSD17B10 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_004493](#).

Gene Aliases

ABAD, ERAB, HCD2, MHBD, HADH2, SCHAD, 17b-HSD10

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

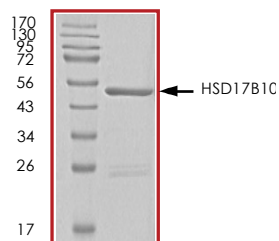
Scientific Background

HSD17B10 gene encodes the protein 17-beta-hydroxysteroid dehydrogenase 10 that is a member of the short-chain dehydrogenase/reductase superfamily SCHAD (1). HSD17B10 gene product is a mitochondrial protein that is involved in lipid metabolism, fatty acid oxidation and steroid hormone metabolism. HSD17B10 protein has been implicated in the development of Alzheimer's disease and mutations in the gene are the cause of 2-methyl-3-hydroxybutyryl-CoA dehydrogenase deficiency (MHBD). Furthermore, HSD17B10 may act as a direct molecular link between beta-amyloid and mitochondrial toxicity (2).

References

1. Yang, S.-Y. et al: Multiple functions of type 10 17-beta-hydroxysteroid dehydrogenase. Trends Endocr. Metab. 16: 167-175, 2005.
2. Lustbader, J. W. et al: ABAD directly links A-beta to mitochondrial toxicity in Alzheimer's disease. Science 304: 448-452, 2004.

Purity



The purity of HSD17B10 was determined to be **>95%** by densitometry. Approx. MW **51kDa**.

HSD17B10 Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog Number	H23-30G
Specific Lot Number	U316-2
Purity	>95%
Concentration	0.2 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.