

TDO2 Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog # T213-30G

Lot # A1351-3

Product Description

Full-length recombinant human TDO2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [NM_005651](#).

Gene Aliases

TDO; TO; TPH2; TRPO

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM 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.

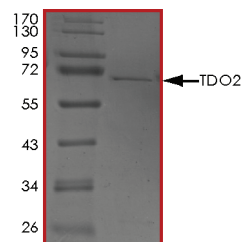
Scientific Background

TDO2 or tryptophan 2, 3-dioxygenase is a heme enzyme that plays a critical role in tryptophan metabolism by catalyzing the first and rate-limiting step of the kynurenine pathway (1). TDO2 may play a role in cancer and immune biology through the suppression of antitumor immune responses, and single nucleotide polymorphisms in TDO2 may be associated with autism (2).

References

1. Comings, D. E. et.al: Localization of human tryptophan oxygenase to 4q31: possible relevance to alcoholism, depression and Tourette syndrome. (Abstract) Am. J. Hum. Genet. 45 (suppl.): A135 only, 1989.
2. Opitz, C. A.et.al: An endogenous tumour-promoting ligand of the human aryl hydrocarbon receptor. Nature 478: 197-203,2011.

Purity



The purity of TDO2 protein was determined to be **>95%** by densitometry. Approx. MW **69 kDa**.

TDO2 Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog #	T213-30G
Lot #	A1351-3
Purity	>95%
Concentration	0.1 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.