

## IDO1 Protein

Full-length recombinant human protein expressed in Sf9 cells

**Catalog # I211-30G**

Lot # A1339-4

### Product Description

Full-length recombinant human IDO1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [BC027882](#).

### Gene Aliases

CD107B, IDO, INDO

### 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.

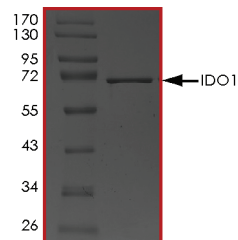
### Scientific Background

IDO1 encodes indoleamine 2,3-dioxygenase (IDO) - a heme enzyme that catalyzes the first and rate-limiting step in tryptophan catabolism to N-formyl-kynurenine which acts on multiple tryptophan substrates including D-tryptophan, L-tryptophan, 5-hydroxy-tryptophan, tryptamine, and serotonin. IDO1 play a main role in a variety of pathophysiological processes such as antimicrobial and antitumor defense, neuropathology, immunoregulation, and antioxidant activity. IDO1 activity in tumor cells may act to impair antitumor responses (1). The efficacy of therapeutic vaccination of cancer patients might be improved by concomitant administration of an IDO inhibitor (2).

### References

1. Logan, G. J. et.al: HeLa cells cocultured with peripheral blood lymphocytes acquire an immuno-inhibitory phenotype through up-regulation of indoleamine 2,3-dioxygenase activity. *Immunology* 105: 478-487, 2002.
2. Uyttenhove, C.et.al: Evidence for a tumoral immune resistance mechanism based on tryptophan degradation by indoleamine 2,3-dioxygenase. *Nature Med.* 9: 1269-1274, 2003.

### Purity



The purity of IDO1 protein was determined to be **>95%** by densitometry. Approx. MW **68 kDa**.

## IDO1 Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog #	I211-30G
Lot #	A1339-4
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.

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

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