

Data Sheet

Mouse IDO1, His-tag

Human, recombinant, N-terminal His-tag

Catalog #: 71196

Lot #: 150521

Conc.: 3.70 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted protein is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. Lob, S. *et al.*, *Cancer Immunol. Immunother.* 2009; **58(1)**: 153-157.
2. Liu, X. *et al.*, *Blood.* 2010; **115(17)**: 3520-3530.
3. Flick, H.E., *et al.*, *Int. Nat. J. Tryptophan Res.* 2013; **6**: 35-45.

Description:

Mouse IDO1, also known as Indoleamine 2,3-dioxygenase 1, GenBank Accession No. NM_008324, a.a. 2-407(end) with an N-terminal His-tag, expressed in the presence of hemin in an *E. coli* cell expression system. MW = 46 kDa.

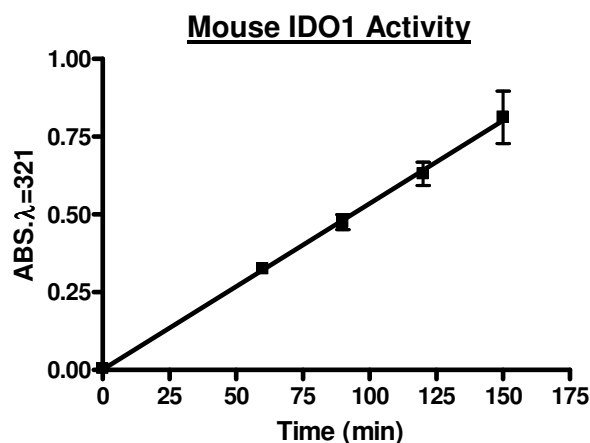
Activity:

Assay Conditions: Assay was performed in 20 mM ascorbate, 10 mM methylene blue, 10 $\mu\text{g/ml}$ catalase, 100 μM NaPi, pH 6.5 containing varying concentrations of tryptophan. 300 ng of IDO1 was added to initiate reaction and aliquots were collected at various time points for Kyn analysis.

Application:

Useful for studying enzyme kinetics, substrate specificity, and screening inhibitors.

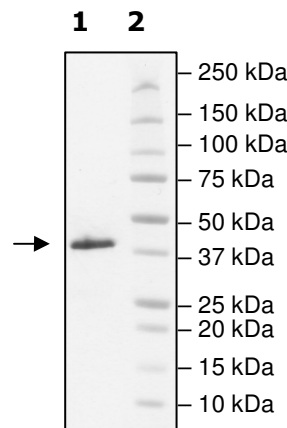
Quality Assurance



**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 2 μg Mouse IDO1
Lane 2:
 Protein Marker

Purity: $\geq 95\%$
MW: 46 kDa



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com