

Data Sheet

IDO1, His-Avi-Tag, Biotin Labeled

Human, Recombinant, N-terminal His-Avi-tag,
Biotin-labeled

Catalog #: 79081

Lot #: 170701-1

Conc.: 1.12 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 0.04% Tween
20, 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. of Tryptophan Res.* 2013; **6**: 35-45.

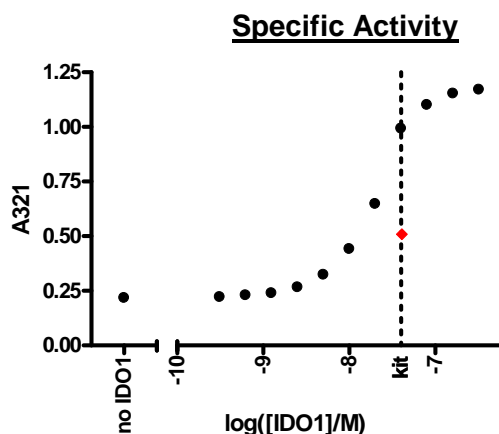
Description:

Human IDO1, also known as Indoleamine 2,3-dioxygenase 1, GenBank Accession No. NM_002164, a.a. 2-403(end) with an N-terminal His-Avi-Tag, expressed in an *E. coli* cell expression system. MW = 48 kDa.

Assay Conditions: The assay was performed by UV absorption using a recombinant IDO1 or TDO2 and L-Tryptophan substrate. The UV absorption signal at 321 nm is correlated with the amount of N-formylkynurenine reaction product of reaction. All of the reactions were conducted at room temperature. The 100 μl reaction mixture in IDO or TDO Assay Buffer contains ~ 40 nM IDO1 (200 ng/well for 71182 or 210 ng/well for 79081) or varying concentrations (0-320 nM) of IDO1, 900 tryptophan, and the coupled reaction components. The reaction mixture incubated for 165 min respectively while reading the UV absorption signal. For the negative control (blank), assay buffer was added instead of the enzyme. Absorption signals were measured using a Tecan Infinite M1000 plate reader.

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

Quality Assurance



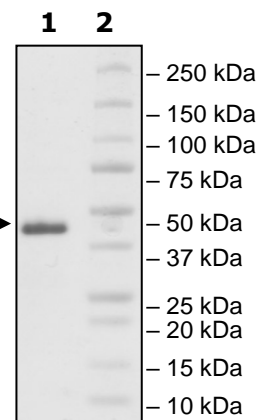
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
IDO1

Lane 2:
Protein Marker

Purity: $\geq 90\%$

MW: 48 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