

PRODUCT INFORMATION



IDO1 (human, recombinant)

Item No. 32559

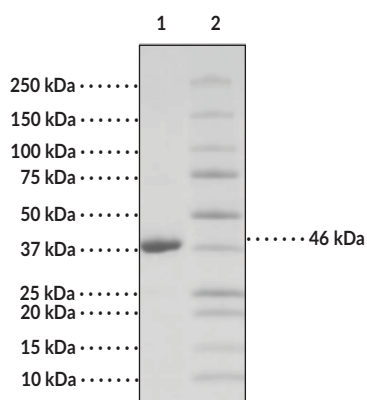
Overview and Properties

Synonyms: Indoleamine 2,3-dioxygenase 1, Indoleamine-pyrrole 2,3-dioxygenase
Source: Active recombinant human N-terminal His-tagged IDO1 expressed in *E. coli*
Amino Acids: 2-403 (full length)
Uniprot No.: P14902
Molecular Weight: 46 kDa
Storage: -80°C (as supplied)
Stability: ≥6 months
Purity: **batch specific** (≥90% estimated by SDS-PAGE)
Supplied in: 40 mM Tris-HCl, pH 8.0, with 110 mM sodium chloride, 2.2 mM potassium chloride, and 20% glycerol

Protein
Concentration: **batch specific** mg/ml
Activity: IDO1 activity is measured by incubating with L-Trp substrate. After a three hour room temperature incubation, activity is determined by measuring the absorption of reaction product at $\lambda = 320-325$ nm

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

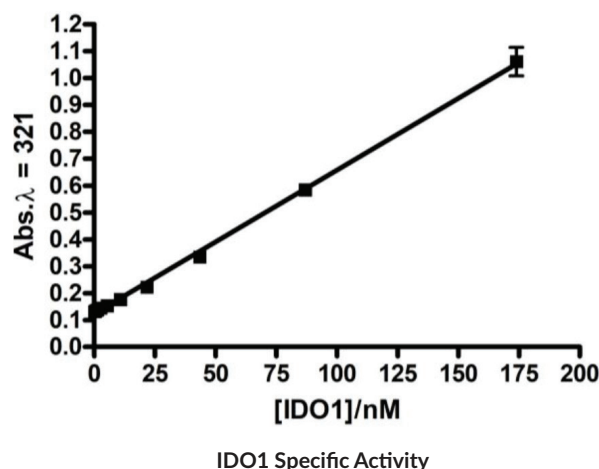
Images



Lane 1: IDO1 (2 µg)
Lane 2: MW Markers

SDS-PAGE Analysis of IDO1.

Representative gel image shown; actual purity may vary between each batch.



WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/10/2020

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Indoleamine 2,3-dioxygenase 1 (IDO1) is a cytosolic heme-containing enzyme involved in tryptophan catabolism that mediates multiple immunomodulatory processes.^{1,2} It exists as a monomer and is composed of a large catalytic domain, which contains the heme group, and a small non-catalytic domain with two immunoreceptor tyrosine-based inhibition motifs (ITIMs) that regulate intracellular signaling.^{3,4} IDO1 catalyzes the oxidation of L-tryptophan (Item No. 29600) to N-formylkynurenine, the first and rate-limiting step of the kynurenine pathway, leading to the production of several metabolites of kynurenine (Item No. 11305), including 3-hydroxy anthranilic acid (Item No. 20512), that have immunoprotective and immunosuppressive effects.¹ IDO1 is widely expressed in several lymphoid and non-lymphoid organs, including lymph nodes, spleen, tonsils, lungs, and intestine, as well as immune cells, including macrophages and dendritic cells (DCs), and endothelial cells.¹ IDO1 activity is regulated at the transcriptional level where expression of IDO1 is increased by the transcription factors NF- κ B, aryl hydrocarbon receptor (AhR), and C/EBP β , which are induced by stimulation with toll-like receptor (TLR) ligands or cytokines and inhibited by hypoxia.¹ IDO1 activity can also be inhibited at the protein level by peroxynitrite (Item No. 81565), leading to IDO1 inactivation, or suppressor of cytokine signaling 3 (SOCS3), which binds to and targets IDO1 for proteasomal degradation. Increased IDO1 activity is associated with numerous pathological conditions, including atherosclerosis, autoimmunity, HIV, and cancer. Cayman's IDO1 (human, recombinant) protein can be used for enzyme assay applications. This protein consists of 402 amino acids and has a calculated molecular weight of 46 kDa.

References

1. Hornyák, L., Dobos, N., Koncz, G., *et al.* The role of indoleamine-2,3-dioxygenase in cancer development, diagnostics, and therapy. *Front. Immunol.* **9**, 151 (2018).
2. Röhrig, U.F., Majjigapu, S.R., Vogel, P., *et al.* Challenges in the discovery of indoleamine 2,3-dioxygenase 1 (IDO1) inhibitors. *J. Med. Chem.* **58**(24), 9421-9437 (2015).
3. Coletti, A., Greco, F.A., Dolciemi, D., *et al.* Advances in indoleamine 2,3-dioxygenase 1 medicinal chemistry. *Med. Chem. Commun.* **8**(7), 1378-1392 (2017).
4. Cheong, J.E. and Sun, L. Targeting the IDO1/TDO2-KYN-AhR pathway for cancer immunotherapy – challenges and opportunities. *Trends Pharmacol. Sci.* **39**(3), 307-325 (2018).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM