

PRODUCT INFORMATION



ARG1 (human, recombinant)

Item No. 32558

Overview and Properties

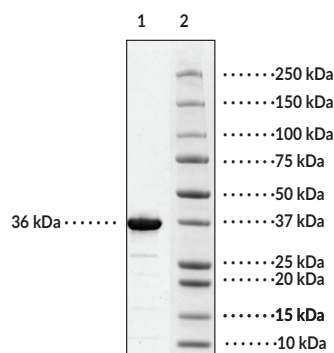
Synonyms: Arginase1, Liver-type Arginase, Type I Arginase
Source: Active recombinant human C-terminal His-tagged ARG1 expressed in HEK293 cells
Amino Acids: 1-322 (full length)
Uniprot No.: P05089
Molecular Weight: 36 kDa
Storage: -80°C (as supplied)
Stability: ≥6 months
Purity: *batch specific* (≥90% estimated by SDS-PAGE)
Supplied in: 8 mM Phosphate buffer, pH 7.4, with 110 mM sodium chloride, 2.2 mM potassium chloride, and 20% glycerol

Protein

Concentration: *batch specific* mg/ml

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

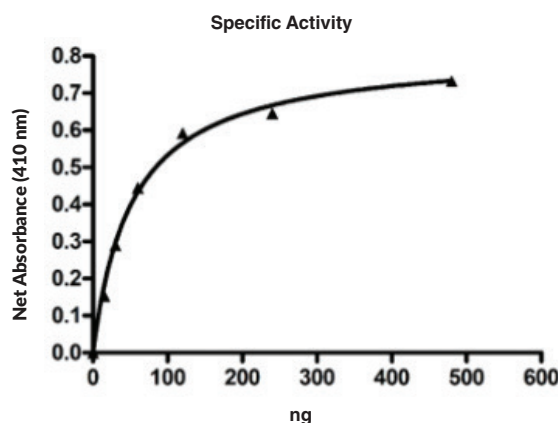
Images



Lane 1: ARG1 (4 µg)
Lane 2: MW Markers

SDS-PAGE Analysis of ARG1.

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, 04/29/2025

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

Arginase-1 (ARG1) is a manganese-metalloenzyme that catalyzes the irreversible hydrolysis of L-arginine into urea and L-ornithine with roles in polyamine synthesis and tissue repair.¹⁻³ It functions as a homotrimer with each subunit containing a binuclear manganese cluster at its active site.² ARG1 is expressed primarily in the liver, and at lower levels in various extrahepatic tissues, including the heart, lung, intestine, and stomach, as well as the CNS, blood vessels, and M2 macrophages.¹ Loss-of-function mutations in ARG1 are associated with the autosomal recessive disorder arginase-1 deficiency that is characterized by neurological impairment, growth retardation, spastic paraparesis, and hyperargininemia.² ARG1 expression is increased in lung homogenates isolated from patients with asthma.⁴ Cayman's ARG1 (human, recombinant) protein can be used for enzyme assay applications.

References

1. Moretto, J., Girard, C., and Demougeot, C. The role of arginase in aging: A systematic review. *Exp. Gerontol.* **116**, 54-73 (2019).
2. Sin, Y.Y., Baron, G., Schulze, A., et al. Arginase-1 deficiency. *J. Mol. Med. (Berl.)* **93(12)**, 1287-1296 (2015).
3. Greenhalgh, A.D., Passos Dos Santos, R., Zarruk, J.G., et al. Arginase-1 is expressed exclusively by infiltrating myeloid cells in CNS injury and disease. *Brain Behav. Immun.* **56**, 61-67 (2016).
4. North, M.L., Khanna, N., Marsden, P.A., et al. Functionally important role for arginase 1 in the airway hyperresponsiveness of asthma. *Am. J. Physiol. Lung Cell Mol. Physiol.* **296(6)**, L911-L920 (2009).

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