

Arginase I (human), (recombinant)

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|--------------|------|
| ALX-201-0810 | 20µg |
|--------------|------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name

ARG1

Formulation

Liquid. In 10mM TRIS-HCl, pH 7.5, containing 1mM β -mercaptoethanol, 1mM $MnCl_2$ and 50% glycerol.

Purity

$\geq 90\%$ (SDS-PAGE)

Source

Produced in *E. coli*. Full length human arginase I.

Specific Activity

$\sim 2.0U/\mu g$ protein. One unit is defined as the amount of enzyme that converts $1\mu mol$ of L-arginine to L-ornithine and urea per min. at 37°C, pH 9.5 (R.T. Schminke, et al.; J. Biol. Chem. **238**, 1012 (1962)).

UniProt ID

P05089

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)