

Bone morphogenetic protein 2 (human), (recombinant)

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-205-C010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	The reconstituted solution can be further diluted in sodium acetate solution and stored at +4°C for up to two weeks. Below 10µg/ml, 0.5% BSA or serum should be added.
Handling	After reconstitution, prepare aliquots and store at -80°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	BMP-2
Biological Activity	Induces alkaline phosphatase expression at an EC ₅₀ = 14nM in C2C12 cells.
Concentration	100-500µg/ml after reconstitution.
Endotoxin Content	<0.02EU/mg protein.
Formulation	Lyophilized.
MW	~25.9kDa.
Purity	≥98% (SDS-PAGE, UV, MS)
Reconstitution	Reconstitute with 20-100µl 50mM sodium acetate solution, pH 4.3.
Source	Produced in <i>E. coli</i> and renatured from inclusion body material.
UniProt ID	P12643

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-be@enzolifesciences.com)

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