

BMPR-1A (human), (recombinant)

Bone morphogenetic protein receptor type-1A (BMPR-1A) is a Type I receptor that complexes with BMP Type II receptors to activate SMAD signaling. BMPR-1A is a receptor for BMP-2, BMP-4, GDF-5, and GDF-6. This product contains the first 152 amino acids of the extracellular portion of BMPR-1A and can be used to interfere with or neutralize ligand signaling.

Ordering Information

[Order Online »](#)

ENZ-PRT260-0020	20µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

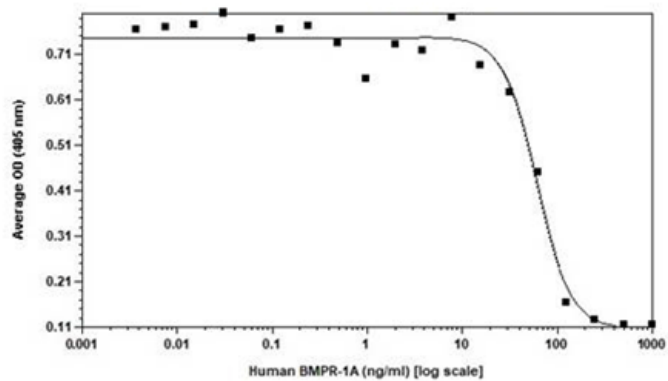


Figure 1: Human BMPR1A inhibition of BMP-4 activity on ATDC5 cells

Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C. For long term storage, it is recommended to dilute to working aliquots in a 0.1% BSA solution.
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	BMP type-1A receptor, Activin receptor-like kinase 3, SKR5
Endotoxin Content	≤1 EU/μg
Formulation	Lyophilized from a sterile (0.2 micron) filtered aqueous solution containing PBS.
MW	Dimer, 40.4/80.8 kDa (362/724 amino acids)
Purity	>95% (SDS-PAGE)
Reconstitution	Reconstitute in sterile water at 0.1 mg/mL.
Source	Produced in genetically modified CHO cells. Contains a C-terminus linker (IEGRMD) and an Fc tag (Human IgG1).
UniProt ID	P36894

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)