

Retinoid X receptor β (human), (recombinant)

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE127-0050	50 μ g
----------------	------------

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	RXRβ
Concentration	2 mg/ml total protein (~0.6 mg/ml RXRβ)
Formulation	Liquid. In 20 mM HEPES, pH 7.9, containing 0.1 M KCl, 0.2 mM EDTA, 0.5 mM DTT and 20% glycerol.
MW	51 kDa
Purity Detail	Partially purified.
Source	Produced in Sf9 insect cells. Full length RXRβ (rH2RIIBP). Produced in a baculovirus expression system.
UniProt ID	P28702

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)