

REP-1 His

Ordering Information

[Order Online »](#)

JBS-PR-105

50µg

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 Rab Escort Protein 1

Formulation liquid (Supplied in 25 mM HEPES pH 7.2, 40 mM NaCl and 5 mM DTT)

MW 75 kDa

Purity > 90 % (SDS-PAGE)

Source fish, recombinant, E. coli

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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