

RaIGDS-RBD^{GST}

Ordering Information

[Order Online »](#)

JBS-PR-365	250µg
------------	-------

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 Ras Binding Domain of the Ral Guanine Nucleotide Dissociation Stimulator

Formulation liquid (Supplied in 30 mM Tris-HCl pH 7.5, 100 mM NaCl, 5 mM MgCl₂ and 5 mM DTE)

MW 40 kDa

Purity > 90 % (SDS-PAGE)

Source human, recombinant, E. coli, Recognized GTPases: Rap-1A, Rap-1B, Rap-2A, Rap-2B, Rit, Rin

Specificity Recognizes GTPases: Rap-1A, Rap-1B, Rap-2A, Rap-2B, Rit, Rin.

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](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)