

[pTyr1158/1162/1163]Insulin receptor and [pTyr1131/1135/1136]IGF- 1 Receptor polyclonal antibody

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

[Order Online »](#)

BML-SA392-0020	20µl
BML-SA392-0100	100µl

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 IR and IGF-1R (Insulin-like growth factor-1 receptor)

Application WB

Formulation Liquid. In Dulbeccoo's PBS, pH 7.3, with 1mg/ml BSA, 50% glycerol and 0.05% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to a portion of human insulin receptor phosphorylated at Tyr1158, Tyr1162 and Tyr1163.

Species Reactivity Human, Mouse

Specificity Recognizes insulin receptor/IGF-1R phosphorylated at Tyr^{1158/1162/1163}.

UniProt ID P06213

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



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)