

Insulin receptor (human), (recombinant) (His- tag) (GST-tag)

INSR (Insulin receptor) (CD220), is a transmembrane receptor activated by insulin. INSR exists as a tetramer consisting of two alpha and two beta subunits linked by disulfide bonds. INSR associates with downstream mediators upon binding to insulin, including IRS1 (insulin receptor substrate 1) and phosphatidylinositol 3'-kinase (PI3K). INSR is expressed in the peripheral nerve, fibroblasts, kidney, liver, skin and striated muscle. Defects in INSR are associated with insulin resistance, Donohue syndrome, noninsulin-dependent diabetes mellitus, and Rabson-Mendenhall syndrome.

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. Prepare aliquots for optimal storage.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CD220, HHF5, INSR
Application Notes	Useful for enzyme kinetics and inhibition studies and for phosphorylation of target substrates.
Formulation	Liquid. In 50 mM Tris, pH 7.4, containing 100 mM NaCl, 0.3 mM DTT and 20% glycerol.
MW	72.3 KDa
Purity	≥90% (SDS-PAGE)
Source	Produced in insect cells using Baculovirus expression system. The cytoplasmic domain of the human insulin receptor (aa 989-1382) is fused with an N-terminal polyhistidine-tagged GST tag at the N-terminus.
UniProt ID	P06213

Last modified: May 29, 2024



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