

# 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.

## Ordering Information

[Order Online »](#)

ENZ-PRT207-0020	20µg
-----------------	------

## 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



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-  
fr@enzolifesciences.com](mailto:info-fr@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)