[pTyr1158/1162/1163]Insureceptor and [pTyr1131/1135/1136]IGF-1 Receptor polyclonal

antibolis Gyed by our Worry-Free Guarantee.

Ordering Information

Order Online »

BML-SA392-0020	20μΙ
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

.

Last modified: May 29, 2024

