Nerve growth factor receptor (human):Fc (human), (recombinant)

Ordering Information

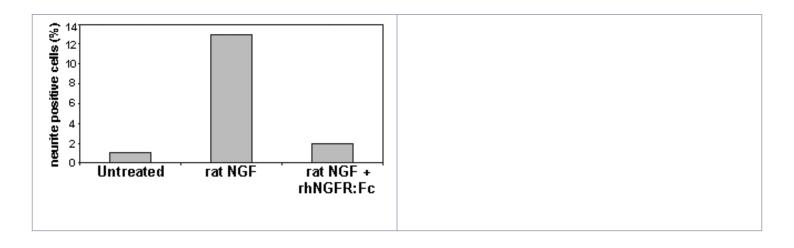
Order Online »

ALX-522-044-C050

50µg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability Stable for at least 6 months after receipt when stored at -20°C. After reconstitution, it is

recommended to make appropriate aliquots and to store them at -20°C.

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name NGFR, TNFRSF 16, Low-affinity nerve growth factor

receptor, p75NTR, p75 neurotrophin receptor, CD271

Concentration 1mg/ml after reconstitution.

Endotoxin Content <0.1EU/μg purified protein (LAL test; Bio Whittaker).

Formulation Lyophilized. Contains PBS.

MW ~80kDa (under reducing conditions).

Purity ≥90% (SDS-PAGE)

Reconstitution Reconstitute in 50µl sterile water. Further dilutions can be

made with medium containing 5% fetal calf serum.

Source Produced in HEK 293 cells. The cysteine-rich domain of

human NGFR (nerve growth factor receptor; p75 NTR) (aa

1-250) is fused to the Fc portion of human IgG1.

SpecificityBinds human, mouse and rat NGF (nerve growth factor).

Technical Info / Product Notes Historical data has shown that NGFR inhibits rat NGF-

mediated activation of rat PC12 cells.

UniProt ID P08138

Last modified: May 29, 2024



info-

eu@enzolifesciences.com