GITR (mouse):Fc (human), (recombinant)

Citations: 5

View Online »

Ordering Information

Order Online »

ALX-522-024-C050

50µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability Stable for at least 6 months after receipt when stored at -20°C.

Handling Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name AITR, TNFRSF 18, Glucocorticoid-induced TNFR-related

protein, Activation-inducible TNFR family receptor

Concentration 1mg/ml after reconstitution.

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

Formulation Lyophilized. Contains PBS.

MW ~45kDa (SDS-PAGE).

Purity ≥95% (SDS-PAGE)

Reconstitution Reconstitute with 50µl sterile water. Further dilutions

should be made with medium containing 5% fetal calf

serum or a carrier protein.

Source Produced in HEK 293 cells. The cysteine-rich region

of mouse GITR (aa 20-152) is fused to the Fc portion of

human IgG1.

Specificity Binds mouse GITRL.

UniProt ID O35714

