

Data Sheet

GITRL

Human, recombinant, N-terminal His-tag

Catalog #: 71190

Lot #: 140826-D

Conc.: 1.43 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol

Stability: >6 months at -80°C . Avoid freeze/thaw cycles. Protein may be diluted to $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .

Endotoxin Level: <1 EU/ μg

References:

1. Shevach, E.M. *et al.*, *Nat. Rev. Immunol.* 2006; **6**: 613-618.
2. Hanabuchi, S. *et al.*, *Blood.* 2006; **107(9)**: 3617-3623.

Description:

Human glucocorticoid-induced TNF-related ligand (GITRL), also known as AITRL, CD357L, or tumor necrosis factor ligand superfamily member 18 (TNFSF18). GenBank Accession No. NM_005092, a.a. 72-199 with an N-terminal His-tag, expressed in a HEK293 cell expression system. MW = 17 kDa. Note: this protein runs at a higher MW by SDS-PAGE due to glycosylation.

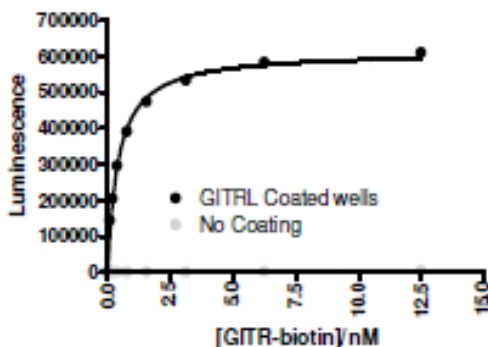
Assay Description: GITR-biotin binding to GITRL was confirmed using the GITRL: GITR Biotinylated] Inhibitor Screening Assay Kit (BPS Cat. #72061).

Application:

Useful for studying protein binding and for screening small molecules and antibodies.

Quality Assurance

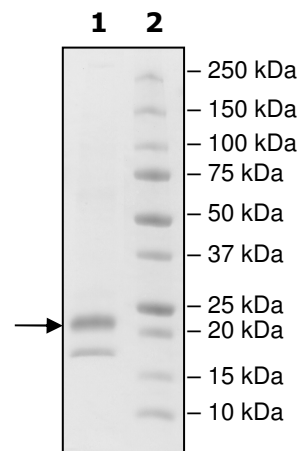
GITRL: GITR-biotin Binding



**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2 μg GITRL
Lane 2:
Protein Marker

MW: 17 kDa + glycans
Purity: $\geq 90\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com