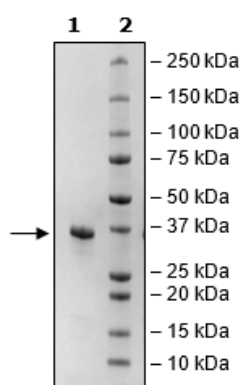


Product Information

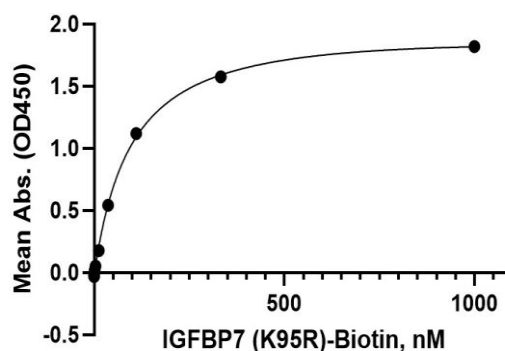
| | |
|---------------------------|--|
| Construct: | IGFBP7 (K95R) (27-282(end)-Avi-His)-(Biotin) |
| Mutation: | K95R |
| Label: | This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%. |
| Concentration: | 5.61 mg/ml |
| Species: | Human |
| Formulated In: | 8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol |
| Expression System: | HEK293 |
| Format: | Aqueous buffer solution |
| Stability: | At least 6 months at -80°C. Avoid freeze/thaw cycles. |
| Storage: | -80°C |
| Genbank Accession: | NM_001553.3 |
| MW: | 30 kDa + glycans |
| Glycosylation: | This protein runs at a higher MW by SDS-PAGE due to glycosylation. |
| Purity: | ≥90% |
| Assay Conditions: | CD93 protein (BPS Bioscience #101543) was coated onto a 96-well plate overnight at 4°C (50 µl/well at a concentration of 4 µg/ml in PBS). The plate was washed 3 times with Immuno Buffer 1 (BPS Bioscience #79311) and blocked using 100 µl of Blocking Buffer 2 (BPS Bioscience #79728) for 1 hour at room temperature. After removing the blocking buffer, 50 µl/well of purified biotinylated IGFBP7 (K95R) protein (BPS Bioscience #101548), serially diluted in Blocking Buffer 2, was added for 30 minutes at room temperature. After 3 more washes, the plate was incubated with Streptavidin-HRP, washed, and incubated with the Colorimetric HRP substrate. The reaction was stopped, and absorbance was read at 450 nm. The Blank value was subtracted from all values. |
| Applications: | Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling. |

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

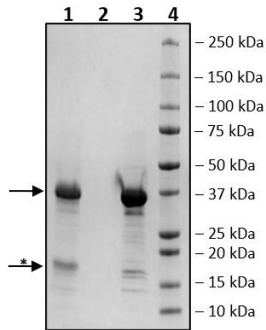


CD93:IGFBP7 (K95R)-Biotin Binding Assay



Quality Control Data

Biotin-Avidin Pulldown



1. Beads
2. Flow thru
3. Control
4. Standards

* Avidin from beads.