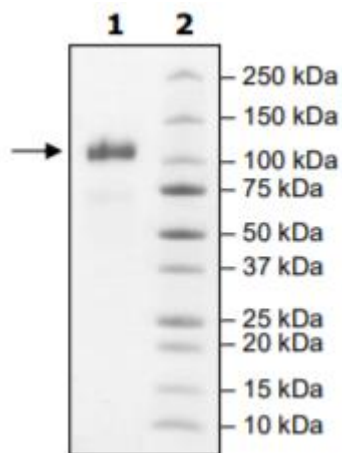


Product Information

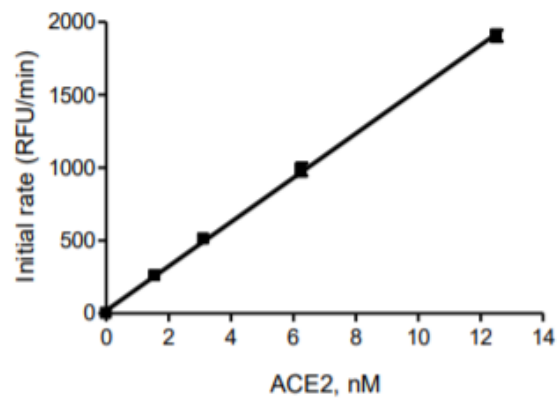
| | |
|---------------------------|--|
| Construct: | ACE2 (18-740-His) |
| Concentration: | 0.14 mg/ml |
| Species: | Human |
| Formulated In: | 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200 mM Imidazole and 20% glycerol. |
| Expression System: | HEK293 |
| Format: | Aqueous buffer solution |
| Stability: | At least 6 months at -80°C. Avoid freeze/thaw cycles. |
| Storage: | -80°C |
| Genbank Accession: | NM_021804 |
| MW: | 85 kDa + glycans |
| Glycosylation: | This protein runs at a higher MW by SDS-PAGE due to glycosylation. |
| Purity: | ≥90% |
| Assay Conditions: | Assay was done according to ACE2 Inhibitor Screening Assay Kit (BPS Bioscience #79923) with ACE2 titrated from 20 ng/rxn to 0.31 ng/rxn. |
| Applications: | As an exopeptidase, ACE2 is useful for the study of enzyme kinetics and to screen for inhibitors. As a cell receptor for coronaviruses infection, including 2019-nCoV, ACE2 is useful for the study of protein-protein interactions. Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling. |

Quality Control Data

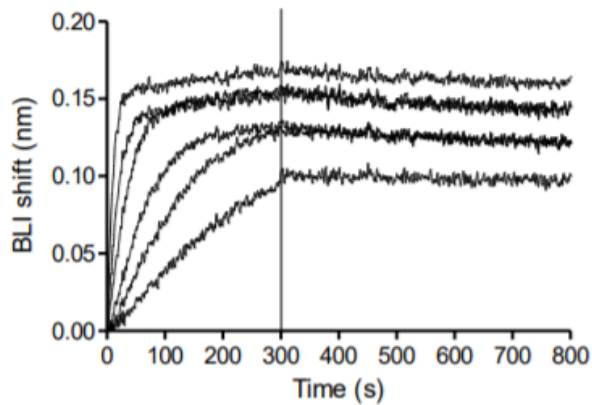
4-20% SDS-Page Coomassie Staining



ACE2 Activity

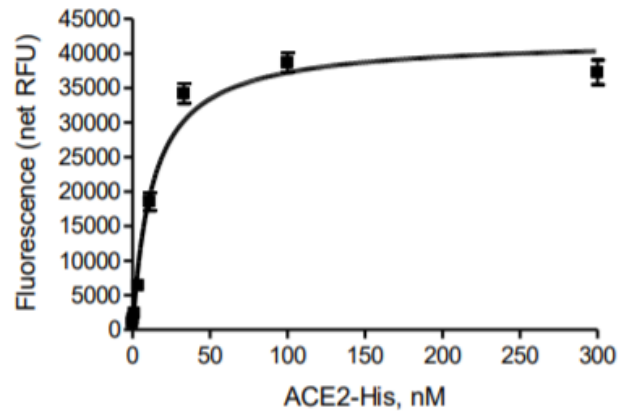


2019-nCoV Spike Protein/Human ACE2, His-tag BLI



Loaded Human ACE2 His-tag (BPS Bioscience #11003) on Anti-His biosensor can bind 2019-nCoV Spike Protein (RBD, mFc Tag) (Sino, 40592-V05H) with an affinity constant of 2 nM as determined in BLI assay (Life Probe Gator).

2019-nCoV Spike Protein/Human ACE2, His-tag DELFIA



Immobilized 2019-nCoV Spike Protein (RBD, mFc Tag) (Sino, 40592-V05H) at 4 $\mu\text{g/ml}$ (200 $\mu\text{l/well}$) can bind human ACE2, His-tag (BPS Bioscience #11003) with a linear range of 1-50 nM. Bound human ACE2, His-tag was quantified with MAb Anti 6HIS-Tb cryptate (Cisbio, 61HISTLB), measuring fluorescence intensity at 340 nm excitation/550 nm emission.