
ACE2, Fc Fusion (Monkey)

Catalog #: 100701

Lot #: 200406

Purity: ≥ 90%

Molecular Weight: 119 kDa

Description: Rhesus monkey angiotensin I converting enzyme 2 (ACE2), also known as ACEH, Genbank Accession No.: ACI04553.1, a.a. 18-739, fused at the C-terminus of the Fc portion of human IgG1, expressed in a HEK293 expression system, MW= 119 kDa. This protein runs at a higher MW due to glycosylation.

Endotoxin Level: <1 EU/μg

Format: Lyophilized from 0.2 μm filtered PBS, pH 7.4

Reconstitution: A quick spin of the vial followed by reconstitution in distilled water to a concentration not less than 0.1 mg/ml. This solution can then be diluted into other buffers.

Storage/Stability: At least 2 years from date of receipt at -20° C.

Quality Control: The purity was determined by SDS-Page and HPLC.

References:

1. Rockx, B., *et al. PLoS One.* **6(4):** e18858 (2011).
2. Li, W., *et al. J. Virology.* **80(9):** 4211-4219 (2006).

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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com