

---

## ACE2, His-tag (Monkey)

**Catalog #:** 100702

**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, expressed in a HEK293 expression system, with a C-terminal His-tag. 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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)