

Data Sheet

CA2

Human, recombinant, C-terminal His-tag

Catalog #: 71094

Lot#: 130509 **Conc.:** 2.85 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 106 mM imidazole and 20% glycerol

Stability: >6 months at -70°C. Avoid freeze/thaw cycles.

References:

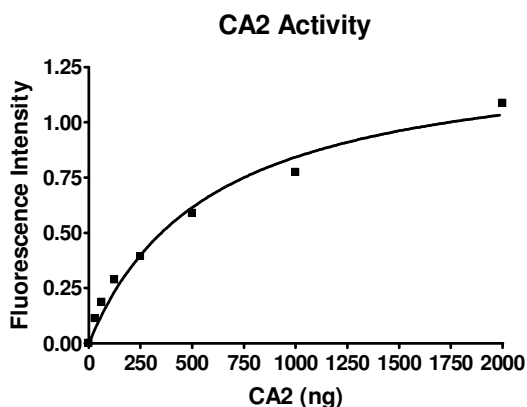
1. Loisel, F.B., *et al.*, *Am J Physiol Cell Physiol.* 2004 Jun; **286(6)**:C1423-33.
2. Shah, G.N., *et al.*, *Hum Mutat.* 2004 Sep; **24(3)**:272.

Description: Human carbonic anhydrase II, also known as CA2, (GenBank Accession No. NM_000067), full length (a.a. 1-260(end)) with C-terminal His-tag, MW = 30 kDa, expressed in an *E. coli* cell expression system.

Specific Activity: 1,400 pmol/min/μg
 Assay conditions: 15 mM Tris-HCl, pH 7.4, 100 mM NaCl, 3 mM p-Nitrophenyl acetate (PNPA) and 0-2 μg CA2. Incubate for 30 minutes at room temperature. Absorbance is measured at 405 nm.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

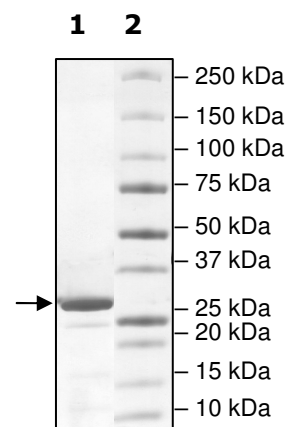
Quality Assurance



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

Lane 1:
 3 μg CA2
Lane 2:
 Protein Marker

MW: 30 kDa
Purity: ≥88%



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



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

**MSHHWGYGKHNGPEHWHKDFPIAKGERQSPVDIDHTAKYDPSLKPLSVSYDQATSL
RILNNGHAFNVEFDDSQDKAVLKGGPLDGTYRLIQFHFHWGSLDGGQSEHTVDKKKYA
AELHLVHWNTKYGDFGKAVQQPDGLAVLGIFLKVGSAPGLQKVVDVLDSIKTKGKSA
DFTNFDPRGLLPESLDYWTYPGSLTTPPILLECVTWIVLKEPISVSSEQVLKFRKLNFNGE
GEPEELMVDNWRPAQPLKNRQIKASFKHHHHHH**

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