

Data Sheet

CA9, His-Tag

Human, Recombinant, C-terminal His-tag

Catalog #: 71101**Lot#:** 130528-2 **Conc.:** 0.21 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 266 mM imidazole and 20% glycerol

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

References:

1. Hilvo, M., *et al.*, *J Biol Chem*. 2008 Oct 10;**283(41)**:27799-809.
2. Chadwick, A.L., *et al.*, *Cell Cycle*. 2013 Jul 30;**12(16)**. [Epub ahead of print]

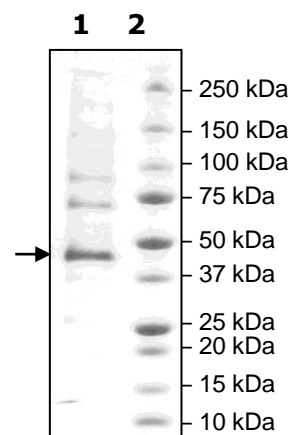
Description: Human carbonic anhydrase IX (CA IX), also known as CA9, (GenBank Accession No. NM_001216), a.a. 59-414 with C-terminal His-tag, MW = 40 kDa, expressed in a baculovirus infected Sf9 cell expression system.

Quality Assurance

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

Lane 1:
3.8 μg CA9
Lane 2:
Protein Marker

Purity: $\geq 65\%$
MW: 40 kDa



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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd, Suite 100

San Diego, CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: support@bpsbioscience.com

Sequence:

MPLGEEDLPSEEDSPREEDPPGEEDLPGEEDLPGEEDLPEVKPKSEEEGSLKLEDLPTV
EAPGDPQEPQNNNAHRDKEGDDQSHWRYGGDPPWPRVSPACAGRFQSPVDIRPQLAAF
CPALRPLELLGFQLPPLPELRLRNNGHSVQLTLPPGLEMALGPGREYRALQLHLHWGAA
GRPGSEHTVEGHRFP AEIHVVHLSTAFARVDEALGRPGGLAVLAAFLEEGPEENSAYEQ
LLSRLEEIAEEGSETQVPGLDISALLPSDFSRYFQYEGSLTTPPCAQQVIWTVFNQTVMLS
AKQLHTLSDTLWGP GDSRLQLNFRATQPLNGRVIEASFPAGVDSSPRAAEPVQLNSCLA
AGDHHHHHH

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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com