

## Data Sheet

### CA2, His-Tag

Human, recombinant, C-terminal His-tag

Catalog #: 71094

Lot#: 130509

Conc.: 2.85 mg/ml

**Description:** Human carbonic anhydrase II, (CA2), GenBank Accession No. NM\_000067, a.a. 1-260(full length), with C-terminal His-tag, expressed in an *E. coli* cell expression system. MW = 30 kDa.

**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:** At least 6 months at  $-70^{\circ}\text{C}$ . 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.

## Quality Assurance

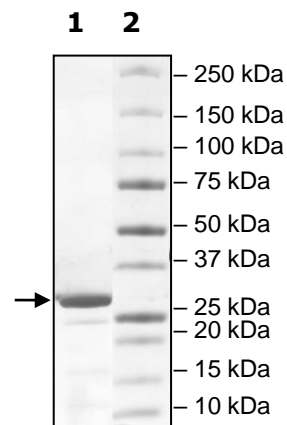
#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
3  $\mu\text{g}$  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 **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

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