

Carbonic anhydrase 2 (human), (recombinant)

Carbonic anhydrase 2 is essential for bone resorption and osteoclast differentiation. It catalyzes the reversible hydration of carbon dioxide. It can hydrate cyanamide to urea. It is also involved in the regulation of fluid secretion into the anterior chamber of the eye. Carbonic anhydrase contributes to intracellular pH regulation in the duodenal upper villous epithelium during proton-coupled peptide absorption. Carbonic anhydrase 2 also stimulates the chloride-bicarbonate exchange activity of SLC26A6

Ordering Information

[Order Online »](#)

ENZ-PRT166-0010	10µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Short Term Storage -20°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CA2, Carbonic anhydrase II, CA-II, Car2

Formulation Lyophilized from 0.22µm filtered solution in 20mM TRIS, pH 8.0, containing 150mM NaCl and 1mM DTT. Normally mannitol or trehalose is added as protectant before lyophilization.

MW ~30 kDa

Purity ≥95% (SDS-PAGE)

Source Produced in HEK 293 cells.

UniProt ID P00918



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)