Cyclooxygenase (sheep), (purified)

Ordering Information

Order Online »

BML-EH001-0001

1mg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name COX

Formulation Liquid. In 0.1M Tris-HCl, pH 7.4 containing 1mM

Tryptophan and 1% Tween-80.

Purity ≥95%

Source Isolated from ram seminal vesicles.

Specific Activity Specific activity with anachidonic acid is 24 units/mg

protein. One unit will consume one micromole of oxygen

per minute at 30°C

UniProt ID P05979 (COX-1), P79208 (COX-2)

Last modified: May 29, 2024



eu@enzolifesciences.com