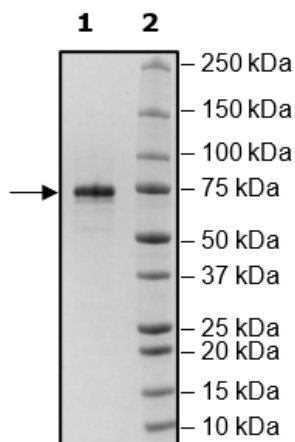


Product Information

Description:	Recombinant human COX1, also known as PTGS1 (prostaglandin-endoperoxide synthase 1), full length encompassing amino acids 1-599 (end). This construct contains a C-terminal FLAG-Tag followed by a His-Tag (6xHis). The recombinant protein was affinity purified.
Species	Human
Construct:	COX1 (1-599(end)-FLAG-His)
Concentration:	1 mg/ml
Expression System:	Sf9
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 0.1% TritonX-100, and 300 μM DDC
MW:	71 kDa
Genbank Accession:	NM_000962
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Assay Conditions:	Activity was measured by titrating various concentrations of enzyme in Assay buffer (100 mM Tris-HCl pH 8, 3 μM EDTA) with 100 μM Ampliflu™ Red and 50 μM arachidonic acid (Cayman Chemical # 90010.1 peroxide free) in a final reaction volume of 100 μL. Fluorescence (ex 535 em 590) was read after 5 minute incubation.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



COX1 Activity

