PARG (bovine thymus)

Citations: 11

Ordering Information

ALX-202-045-UC01

O.1U

Manuals, SDS & CofA

View Online »

Handling & Storage

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

use the following enzyme dilution buffer: 50mM KPO4, pH 7.5, 50mM KCl, 10 mM β-

mercaptoethanol, 10% glycerol, 1mM DTT, 0.1% Triton.

Handling Avoid freeze/thaw cycles.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Poly(ADP-ribose) glycohydrolase

Formulation Liquid. In 50mM K-PO₄ buffer, pH 7.5, containing 50mM

potassium chloride, 0.1% Triton-x 100, 10% glycerol and

1mM DTT, 10mM β-Mercaptoethanol.

Purity Detail Purified by substrate chromatography. The enzyme is

composed of two major bands at 74kDa and 59kDa.

Source Isolated from calf thymus.

Specific Activity ~20'000U/mg protein. 1U is defined as the amount of

enzyme that hydrolyzes 1nmol of poly(ADP-ribose) per

min. at 37°C.

UniProt ID 002776

Last modified: May 29, 2024

