

Product Information

Description:	Fluorogenic PARG Substrate is a fluorogenic ADP-ribose substrate in which the fluorophore is quenched by the presence of the ribose.
Supplied As:	Aqueous solution.
Storage:	Upon receipt, store at -80°C.
Stability:	When stored at -80°C, the product is stable for six months after first thaw. Avoid repeated freeze/thaw cycles.
Application:	It is suitable for use in PARG (poly(ADP-ribose) glycohydrolase) assays using cell extracts or purified proteins. This mixture has been optimized for use with: <ul style="list-style-type: none">• PARG Fluorogenic Assay Kit (BPS Bioscience #78858) Use in other assays may require optimization.

Validation Data

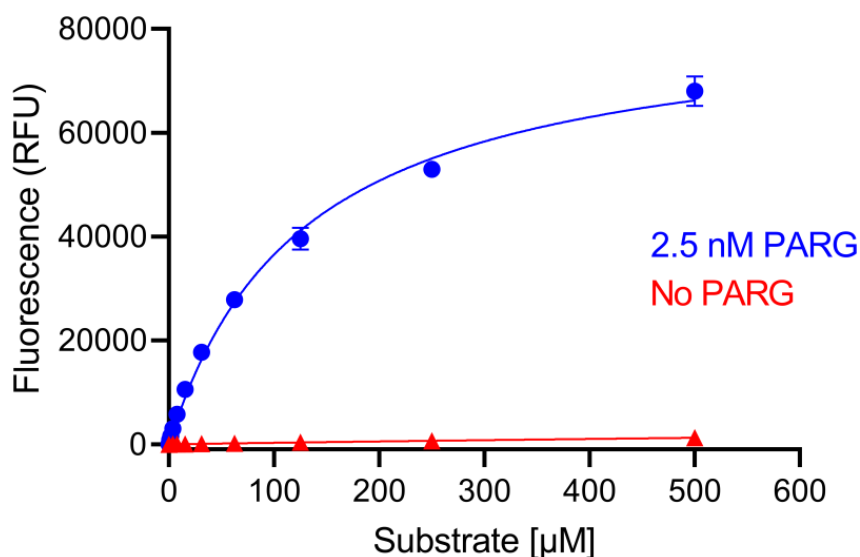


Figure 1: Fluorogenic PARG Substrate allows specific studies on PARG activity.

Fluorogenic PARG Substrate was incubated with no PARG or increasing amounts of purified PARG protein. PARG activity was measured using BPS Bioscience PARG Fluorogenic Assay Kit.