

ADP-ribose-pNP

Catalog number: 11700 Unit size: 1 mg

Component	Storage	Amount
ADP-ribose-pNP	Freeze (<-15 °C), Minimize light exposure	1 mg

OVERVIEW

ADP-ribose-pNP is a colorimetric substrate for assessing activity of poly(ADP-ribose)polymerase (PARP) enzymes. The absorbance of released p-nitrophenol is determined at 405 nm, and the slope of the calibration curve is used to convert the absorbencies to moles of product generated. With ADP-ribose-pNP as the colorimetric substrate, PARP-1 was determined to have the largest Km and Vmax values (151 uM and 1.30 nmolmin-1mg-1 respectively) followed by tankyrase-1 (82 uM and 18 pmolmin-1mg-1 respectively) and VPARP (46 uM and 2 pmolmin-1mg-1 respectively). This colorimetric substrate can be used to determine the kinetic parameters for PARP-1, tankyrase-1, and VPARP, and to screen small-molecule inhibitors of PARP-1, tankyrase-1, and VPARP. ADP-ribose-pNP-based continuous assay has considerable advantages over standard discontinuous PARP assays, enabling the high throughput screening of PARP-1, tankyrase-1, and VPARP activities and their inhibitors.

AT A GLANCE

Important The following recommended procedure can be adapted for measuring PARP-1, tankyrase-1, and VPARP activities and their inhibitors. The optimum conditions must be determined experimentally for each test.

KEY PARAMETERS

Instrument: Absorbance microplate reader

Absorbance: 405 nm
Recommended plate: Solid white

PREPARATION OF STOCK SOLUTIONS

Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20 $^{\circ}$ C after preparation. Avoid repeated freeze-thaw cycles.

1. ADP-ribose-pNP stock solution:

Make 5 - 10 mM stock solution in $\rm H_2O$.

Note The stock solution should be used promptly.

PREPARATION OF WORKING SOLUTION

ADP-ribose-pNP working solution:

Prepare 0.25 mM assay solution by diluting the stock solution with assay buffer $(50 \text{mM Tris}, 10 \text{mM MgCl}_2, \text{pH } 8.0)$.

SAMPLE EXPERIMENTAL PROTOCOL

- Add 0.01 mL/well of sample solution into 0.09 mL/well assay solution to make a final volume of 0.1 mL in a 96-well clear plate.
- 2. Monitor the plate using an absorbance microplate reader at 405 nm.

EXAMPLE DATA ANALYSIS AND FIGURES

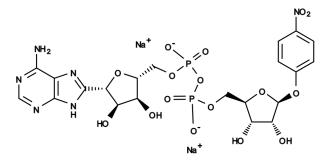


Figure 1. Chemical structure for ADP-ribose-pNP

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.