

AP0002 Protocol

Sandwich ELISA Assay Protocol for Anti-PEG Antibody Pair (AP0002)

Sandwich ELISA Assay Protocol for Anti-PEG Antibody Pair (AP0002):
https://www.novusbio.com/products/polyethylene-glycol-antibody-pair_ap0002
Please centrifuge before opening the vials.

The test protocol is a guideline, user need to determine their optimal experimental condition for best performance.

SECTION 1 Equipments & Reagents

1.1 Anti-PEG Antibody pair

Polyethylene Glycol Matched Antibody Pair (Catalog #: AP0002). This matched antibody pair set binds to the repeating subunits of the polyethylene glycol polymer and can be employed to detect and quantify PEGylated compounds.

1.2 Secondary reagent

Streptavidin-HRP (Jackson ImmunoResearch, Catalog #: 016-030-084)

1.3 Coating buffer (1 Liter)

5.3 g Na₂CO₃ + 4.2 g NaHCO₃, pH=8.0 (adjust pH with 1N NaOH)

1.4 1x PBS

0.14 M NaCl, 2.7 mM KCl, 1.5 mM KH₂PO₄, 8.1 mM Na₂HPO₄, pH 7.4

1.5 Blocking solution

5% skim milk in 1x PBS

1.6 Dilution buffer

2% skim milk in 1x PBS

1.7 PBS-T

1x PBS containing 0.2% Tween-20

1.8 ELISA plates

NUNC MaxiSorp(TM) High Protein-Binding Capacity ELISA plates (Catalog #: 44-2404)

1.9 HRP substrate

50 mg/ml ABTS (Sigma #A-1888) in 100 mM phosphate-citrate buffer pH 4.0 (17.4 g K₂HPO₄, 21 g citric acid in 1 Liter

Q-H₂O). Immediately before use, add 3 ul of 30% H₂O₂ per 10 ml ABTS substrate solution.

SECTION 2 - Assay Protocol

2.1 Dilute capture antibody to 5 ug/ml in coating buffer.

2.2 Add 50 ul diluted capture antibody per well and incubate at 37C for 4 h and then at 4C overnight.

2.3 Wash plates 3 times with 1x PBS.

2.4 Add 200 ul blocking solution per well for 2 hours at room temperature.

2.5 Dilute PEG-compound (analyte) in dilution buffer to suitable concentrations.

2.6 Wash wells 3 times with 1x PBS.

2.7 Add graded concentrations of PEG-compound (50 ul/well) and incubate 2 h at room temperature.

2.8 Wash with PBS-T 3 times and 1x PBS 2 times.

2.9 Add 50 ul/well detection antibody (5 ug/ml in dilution buffer) for 1 h at room temperature.

2.10 Wash wells with PBS-T 3 times and with 1x PBS 2 times.

2.11 Add 50 ul/well streptavidin-HRP (1 ug/ml in dilution buffer), and incubate for 1 h at room temperature.

- 2.12 Wash wells with PBS-T 6 times and with 1x PBS 2 times.
- 2.13 Add 100 ul/well freshly prepared ABTS substrate for 30 min in dark at room temperature.
- 2.14 Read absorbance of the wells at 405 nm.