

## H00002817-A01 Protocol

### Protocol specific for Glypican 1/ GPC1 Antibody (H00002817-A01)

Protocol specific for Glypican 1/ GPC1 Antibody (H00002817-A01): [https://www.novusbio.com/products/glypican-1-antibody\\_h00002817-a01](https://www.novusbio.com/products/glypican-1-antibody_h00002817-a01)  
 ELISA Protocol

1. Coat antigen (200 ng/well) onto the wells in a 96 well microtiter plate.
  2. Block unbound sites with 5% skim milk in PBS.
  3. Apply hybridoma culture supernatant/ ascites/ purified Ig as primary antibody. Incubate the plate at room temperature for two hours.
  4. Wash 4 times with PBST.
  5. Apply HRP conjugated secondary antibody, and incubate the plate at room temperature for one hour.
  6. Wash 8 times with PBST.
  7. Apply 0.1 ml OPD in citric acid buffer, and incubate at room temperature for 20minutes. Read the plate in ELISA reader at 450 nm.
- Between each step, plates were adequately washed using PBST.
  - Secondary antibody dilution, 1:1000.
- Primary Antibody/Dilution:
- Poly sera @ 1500X
  - Cultured Supernatant @ 1X
  - Ascites @ 1000X
  - Purified Ig @ 1 ug/ml
- Diluents:
- 5% skim milk in PBST
- Material:
- PBST, 0.2% Tween 20
  - Citric acid buffer, pH 5.0
  - OPD: Sigma, P-1526

### Western Blot Protocol

1. Antigens were denatured and loaded onto polyacrylamide gel (200 ng/ lane). Run the gel at 150V for 80minutes when samples enter separating gel.
  2. Transfer antigens onto PVDF membrane.
  3. Block PVDF membrane in 5% skim milk in PBST at room temperature for one hour.
  4. Apply hybridoma culture supernatant/ ascites/ purified Ig as primary antibody. Incubate the membrane at room temperature on an orbital shaker for one hour.
  5. Wash 5 times with PBST.
  6. Apply HRP conjugated secondary antibody, and incubate the membrane at room temperature on an orbital shaker for one hour.
  7. Wash 5 times with PBST.
  8. Use UVP autochemi/ ECL system for signal detection (to visualize the result).
- Between each step, plates were adequately washed using PBST.
  - Secondary antibody dilution, 1:10,000.
- Primary Antibody/Dilution:
- Poly sera @ 6000X
  - Cultured Supernatant @ 2X
  - Ascites @ 1000X
  - Purified Ig @ 1ug/ml
- Diluents:
- 5% skim milk in PBST