

## NB600-101 Protocol

### Immunohistochemistry protocol for Calreticulin Antibody (NB600-101)

[[URL:[https://www.novusbio.com/products/calreticulin-antibody\\_nb600-101](https://www.novusbio.com/products/calreticulin-antibody_nb600-101)]][[Caption:Calreticulin Antibody]]  
Immunohistochemistry Procedure

- 1) Paraffin-embedded sections were treated using an HIER [heat-induced epitope retrieval] protocol.
- 2) NB 600-101 was diluted 1:50 with a DAKO antibody diluent
- 3) Sections and primary antibody were incubated at RT for 30 minutes.
- 4) Sections were incubated with a DAKO secondary, as per manufacturer's protocol.
- 5) Sections were then stained with DAB and H&E.
- 6) Sections were mounted with Faramount solution.
- 7) Result of staining was strong.

\*\*NOTE: normal colon mucosa and adenocarcinoma tissues were used as positive controls for this antibody.