



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NB600-101

NB600-101 Protocol

Immunocytochemistry/Immunofluorescence protocol for Calreticulin Antibody (NB600-101)

[[URL:https://www.novusbio.com/products/calreticulin-antibody_nb600-101]][[Caption:Calreticulin Antibody]]
Immunofluorescence Procedure

- 1) Cell cultures were treated using an HIER [heat-induced epitope retrieval] protocol.
- 2) Calreticulin polyclonal antibody [NB 600-101] was diluted 1:50 with a DAKO antibody diluent
- 3) Cultures and primary antibody were incubated at RT for 30 minutes.
- 4) Cultures were incubated with an Alexa green fluorescent secondary, as per manufacturer's protocol.
- 5) Cultures were observed on a Zeiss Axioconvert microscope and with DigiPix software.

**NOTE: COS7 and HCT29 cell cultures were used as positive controls for this antibody.