

Validation data for Anti-hSTING-IgG

<https://www.invivogen.com/anti-hsting>

For research use only

Version 18I28-MM

Anti-hSTING-IgG is a monoclonal mouse IgG1 antibody against human STING (hSTING). It recognizes the most prevalent isoform “wild-type” R232 (R71-G230-**R232**-R293; figure 2D) and HAQ (**H71-A230**-R232-**Q293**; figure 2B) hSTING variants. This antibody is suitable for use in ELISA and Western blot analysis of hSTING.

ELISA

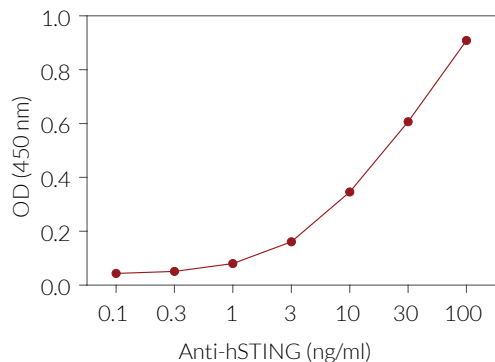


Figure 1. Anti-hSTING-IgG recognition of recombinant hSTING.

Recombinant hSTING was immobilized at 1 µg/mL (50 µL/well) in 96-well plate, followed by the addition of Anti-hSTING capture antibody and HRP-conjugated secondary antibody. Absorbance was read at 450 nm following addition of HRP substrate.

Western blot

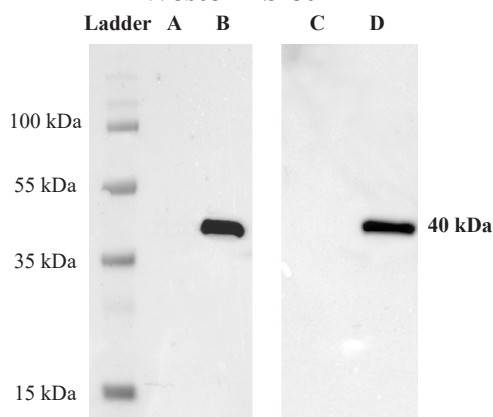


Figure 2. Detection of human STING by Western blot.

Western blot analysis of lysates (40 µg) using Anti-hSTING-IgG (0.2 µg/ml), followed by HRP-conjugated goat anti-mouse IgG1 secondary antibody (1:3000 dilution).

A: THP1-Dual™ KO-STING cells

B: THP1-Dual™ cells (endogenously expresses the HAQ hSTING variant)

C: 293T-Dual™ Null cells (does not express functional hSTING protein)

D: 293T-Dual™ hSTING-R232 cells (stably transfected with the R232 hSTING variant)

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873
InvivoGen USA (International): +1 (858) 457-5873
InvivoGen Europe: +33 (0) 5-62-71-69-39
InvivoGen Hong Kong: +852 3622-3480
E-mail: info@invivogen.com