

Validation data for Anti-hVISTA-hIgG1

<https://www.invivogen.com/anti-hvista-mab>

For research use only

Version 19J24-ED

Anti-hVISTA-hIgG1 is a recombinant monoclonal antibody (mAb) featuring a variable region that recognizes human VISTA and the constant region of the native human IgG1 isotype (hIgG1). VISTA is an important immune checkpoint expressed on myeloid and T cells, that directly suppresses T cell activation and proliferation. The binding of Anti-hVISTA-IgG1 has been validated using engineered VISTA-expressing Raji cells (Figure 1).

Binding of Anti-hVISTA-hIgG1 to target cells

The binding of Anti-hVISTA-hIgG1 to cell-expressed VISTA has been validated using engineered VISTA-expressing Raji cells (Raji-hVISTA) compared to Raji-null cells, which do not express VISTA. A clear shift to the right of the fluorescence peak is noted when Anti-hVISTA-hIgG1 is incubated with Raji-hVISTA (right) and not Raji-null (left) cells, indicating strong and specific binding.

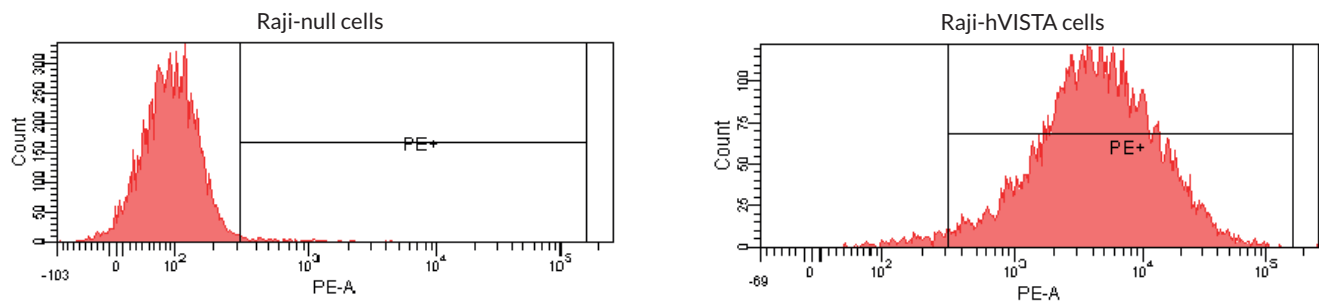


Figure 1: Anti-hVISTA-hIgG1 (1 µg) was added to Raji-null (negative control) and Raji-hVISTA cells (500 000 cells/ml) and incubated at room temperature for 30 minutes. Following this, a secondary antibody, PE, was added and incubated again at room temperature for 30 minutes. The binding affinity was then measured using flow cytometry.

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 3-622-34-80

E-mail: info@invivogen.com