

Validation data for Anti-mCD25-PC-mIgG2a InvivoFit™

<https://www.invivogen.com/anti-mcd25-migg2a-invivofit>

For research use only

Version 21K15-NJ

Anti-mCD25-PC-mIgG2a InvivoFit™ is an anti-mCD25 mAb featuring the variable region of the previously described rat IgG1 clone PC-61.5.3 and a murinized IgG2a constant region. The binding capacity of Anti-mCD25-PC-mIgG2a InvivoFit™ to the murine CD25 (mCD25) antigen has been validated by ELISA, using coated mCD25 protein and a Anti-mIgG-HRP secondary antibody (Figure 1).

Validation of Anti-mCD25-PC-mIgG2a InvivoFit™ by ELISA

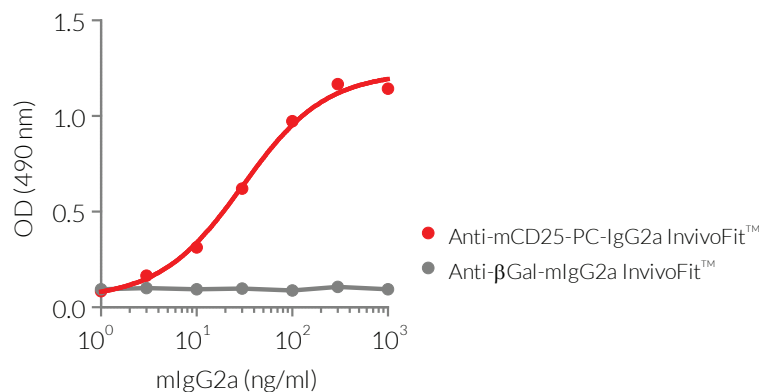


Figure 1: Binding of Anti-mCD25-PC-mIgG2a InvivoFit™ mAb to coated mouse CD25 protein.

Mouse CD25 (2 µg/ml) was coated on ELISA plates overnight. A 3-fold serial dilution of the Anti-mCD25-PC-mIgG2a InvivoFit™ mAb (red curve) or of the Anti-βGal mIgG2a control antibody (grey curve) was performed for the capture step. A HRP-labelled anti-mIgG antibody (1/1000 dilution) and the HRP substrate OPD (o-phenylenediamine dihydrochloride) were used for the detection step. Absorbance was read at 490 nm.

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 Asia: +852 3622-3480
E-mail: info@invivogen.com