## CD162 recombinant monoclonal antibody (PSGL1/8135R)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

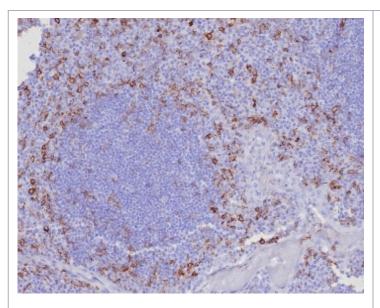
**Order Online** »

**ENZ-ABS877-0100** 

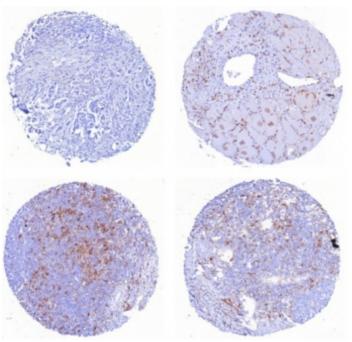
100µg

Manuals, SDS & CofA

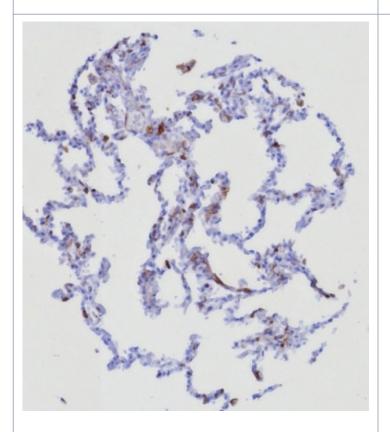
**View Online »** 



IHC staining of FFPE human tonsil tissue with CD162 antibody (clone PSGL1/8135R). HIER: boil tissue sections in pH 9 10mM Tris with 1mM EDTA for 20 min and allow to cool before testing.



IHC staining of FFPE human (1) mammary carcinoma (2) transverse muscle (3) tonsil and (4) tonsil tissue with CD162 antibody (clone PSGL1/8135R) . HIER: boil tissue sections in pH 9 10mM Tris with 1mM EDTA for 20 min and allow to cool before testing.



IHC staining of FFPE human lung tissue with CD162 antibody (clone PSGL1/8135R). HIER: boil tissue sections in pH 9 10mM Tris with 1mM EDTA for 20 min and allow to cool before testing.

## **Handling & Storage**

Handling Aliquot the CD162 antibody and store frozen at -20°C or colder. Avoid freeze/thaw

cycles.

-20°C **Long Term Storage** 

**Shipping** Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

**Alternative Name** P-selectin glycoprotein ligand 1, PSGL-1

**Application** IHC (PS)

Clone PSGL1/8135R

**Formulation** Liquid. 0.2 mg/ml in 1X PBS containing 0.1 mg/ml BSA

(US sourced), 0.05% sodium azide.

Host Rabbit

Recombinant full-length human CD162. **Immunogen** 

Isotype lgGκ

**Purity Detail** Protein A/G affinity purified.

**Recommendation Dilutions/Conditions** Immunohistochemistry (FFPE) (1-2µg/ml for 30 min at

RT)Suggested dilutions/conditions may not be available for

all applications. Optimal conditions must be determined

individually for each application.

**Species Reactivity** Human

**UniProt ID** Q14242

This antibody is covered by our Worry-Free Guarantee **Worry-free Guarantee** 

Last modified: July 12, 2024



info-

eu@enzolifesciences.com