

CD163 (human) clonal antibody (K20-T)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-810-213-R100	100µl
ALX-810-213-R500	500µl

Manuals, SDS & CofA

[View Online »](#)

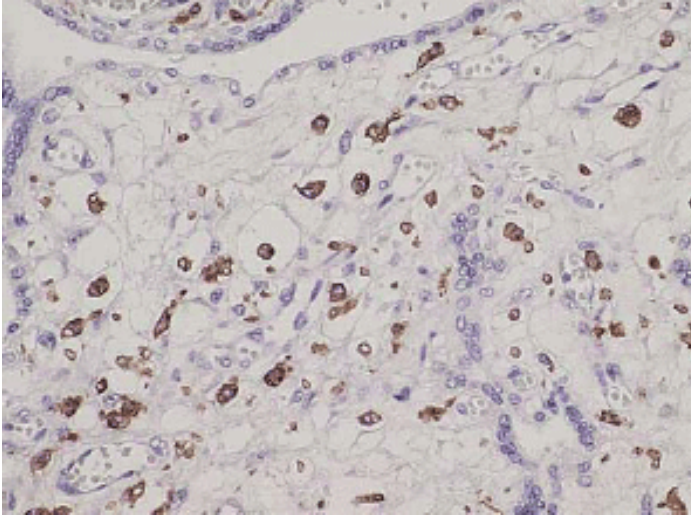


Figure: Formalin-fixed and paraffin-embedded human placenta (4µm) stained with cAb to CD163 (human) (Prod. No. ALX-810-213) shows strong positive cytoplasmic immunostaining of monocytes and macrophages. Kindly performed and provided by Katarína Poliaková, MD and Lubomír Straka, MD, Ph.D. from Clinical Pathology Presov, Ltd., Presov, Slovakia.

Handling & Storage

Use/Stability	Stable for at least 12 months after receipt when stored as recommended.
Handling	After opening, prepare aliquots and store at +4°C. Avoid freeze/thaw cycles. Centrifuge the vial before use.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Scavenger receptor cysteine-rich type 1 protein M130, Hemoglobin scavenger receptor
Application	IHC (FS), IHC (PS)
Clone	K20-T
Formulation	Liquid. In 20mM TRIS/HCl, pH 8.0, containing 20mg/ml BSA and 0.05% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to an N-terminal sequence of human CD163.
Isotype	IgG
Purity Detail	Major clone of IgGs obtained from immunoaffinity purified immunoglobulins corresponding to immunogenic peptide.
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections, 1:100-1:200)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human

Technical Info / Product Notes

Clonal Rabbit Antibody is a pure homogenous fraction of rabbit immunoglobulin (IgG) which corresponds to a unique linear epitope on the target protein. This epitope is designed after a detailed analysis of the antigen structure, regulatory post-translational modifications, protein/protein interaction domain and protein folding. The chosen sequence (peptide) is synthesized and the rabbit is immunized. The clone is then separated from the crude antiserum by “*in vitro* cloning technology”. The resulting product is monospecific and characterized by superior specificity, affinity and avidity.

UniProt ID

Q86VB7

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)