

CD44std (human) monoclonal antibody (SFF-304)

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

Ordering Information

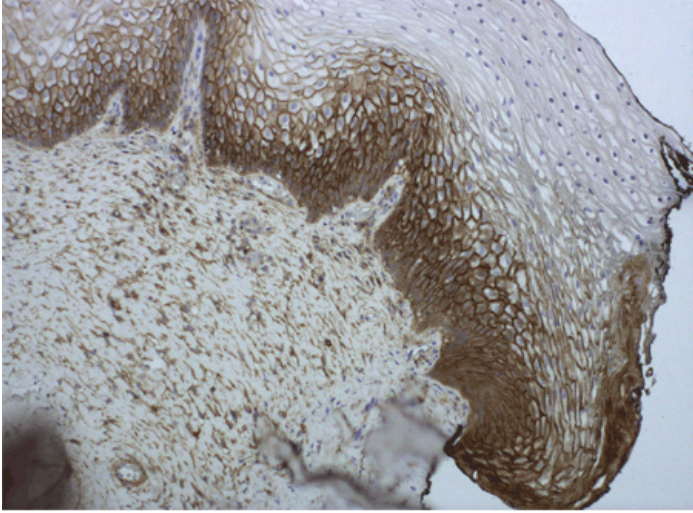
[Order Online »](#)

ALX-801-0890

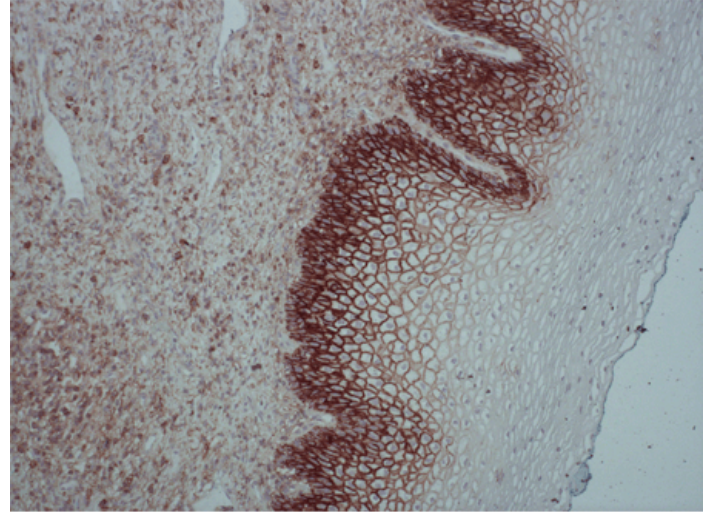
Purified - 100µg

Manuals, SDS & CofA

[View Online »](#)



Cervical tissue assayed with CD44std (human) monoclonal antibody (SFF-304) (ALX-801-089). Developed with POLYVIEW® PLUS HRP (anti-mouse) reagent (ENZ-ACC104) and HIGHDEF® DAB Chromogen/Substrate Set (ENZ-ACC105).



Cervical tissue assayed with CD44std (human) monoclonal antibody (SFF-304) (ALX-801-089). Developed with POLYVIEW® PLUS HRP (anti-mouse) reagent (ENZ-ACC104) and HIGHDEF® DAB Chromogen/Substrate Set (ENZ-ACC105).

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	Flow Cytometry, IHC (FS), IHC (PS), WB
Clone	SFF-304
Formulation	Liquid. In PBS, pH 7.2-7.4.
Host	Mouse
Immunogen	Recombinant human CD44.
Isotype	IgG1
Recommendation Dilutions/Conditions	<p>Flow Cytometry: Recommended concentration for the unlabelled antibody is 2µg/ml as determined on HACAT cells. Immunohistochemistry (frozen sections, paraffin sections): Can be used to stain both indirectly or directly acetone-fixed cryostat sections or cell smears. Recommended is the alkaline phosphatase-anti-alkaline-phosphatase (APAAP) or peroxidase anti-peroxidase (PAP) procedures or the three stage immunoperoxidase technique on acetone-fixed cryostat sections. It may be used at a concentration of 2-10µg/ml for cryostat sections. Suitable for paraffin-embedded tissue after microwave treatment. Western Blot: Recommended concentration: 2-10µg/ml. Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.</p>
Species Reactivity	Human
Specificity	Does not distinguish between CD44 splice variants.

UniProt ID

P16070

Worry-free Guarantee

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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