CD44std (human) monoclonal antibody (SFF-304)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

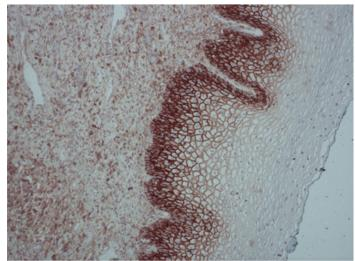
ALX-801-089-C100

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.

lsotype lgG1

Recommendation Dilutions/Conditions 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.

Species Reactivity Human

Specificity Does not distinguish between CD44 splice variants.

P16070

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

