Cytokeratin 8, 18 & 19 monoclonal antibody (2A4)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BML-CA1141-0100

100µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CK8/CK18/CK19

Application ICC, WB

Application NotesDetects bands of 52 kDa, 45 kDa and 40 kDa by Western

Blot.

Clone 2A4

Formulation Liquid. In PBS containing 1% BSA and 0.09% sodium

azide.

Host Mouse

Immunogen Cytokeratins isolated from cultured porcine kidney

epithelial cells (LLC-PK).

lsotype lgG1

Purity Detail Protein A affinity chromatography.

Species Reactivity Human, Mouse, Porcine, Rat

Specificity Recognizes cytokeratin proteins 8, 18 and 19.

UniProt ID P05787 (human CK8), P05783 (human CK18), P08727

(human CK19)

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com