

# 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 Notes Detects 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).

Isotype IgG1

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](#).



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)