

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

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

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-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

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

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

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

Last modified: May 29, 2024