

# Cytokeratin 1 (human) monoclonal antibody (34βB4) (Ready-to- Use)

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

## Ordering Information

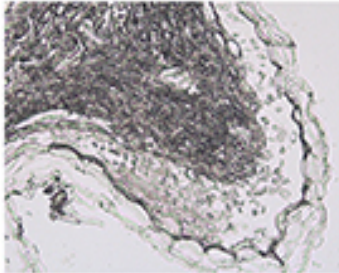
[Order Online »](#)

ENZ-30904	~6 µg - 6ml
-----------	-------------

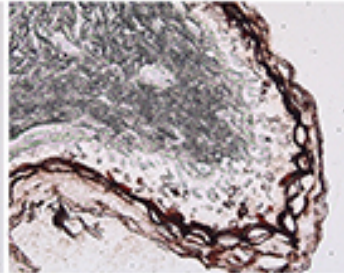
## Manuals, SDS & CofA

[View Online »](#)

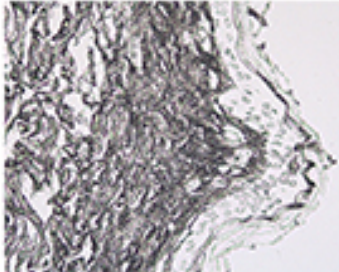
(-) control, 1X Dilution Buffer



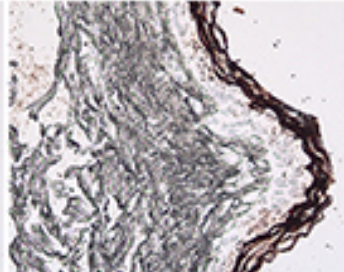
(+) ENZ-C34904



(-) control, 1X Dilution Buffer



(+) ENZ-C34904



Immunohistochemistry analysis of human skin tissue using Cytokeratin 1 (human), mAb (34 $\beta$ B4) at a dilution of 1:1.

# Handling & Storage

Long Term Storage            +4°C

Shipping                        Blue Ice

# Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	CK1
Application	IHC (PS)
Clone	34βB4
Formulation	Liquid. <i>Ready-to-use</i> solution in PBS containing 0.1% sodium azide and 1% BSA.
Host	Mouse
Immunogen	Keratin solubilized from human stratum corneum.
Isotype	IgG1
Purity Detail	Affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections; for optimal results, fixation in alcohol-based fixatives, such as methacarn, is recommended. Pretreatment with proteolytic enzymes such as trypsin may be required when formalin-fixed sections are to be stained. Choose a good quality enzyme-linked anti-mouse IgG1 antibody for detection). Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually.
Source	Purified from ascites.
Species Reactivity	Human

## Specificity

34βB4 is specific for “high molecular weight” cytokeratin 1, 68 kDa, of “complex” epithelium. The antibody reacts with all squamous epithelium and squamous carcinomas. It does not react with adenocarcinomas of either simple epithelium (hepatoma) or ductal origin (breast, pancreas).

## UniProt ID

P04264

## 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](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)