

Cytokeratin 1 (human) monoclonal antibody (34βB4)

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

Citations: 2

[View Online »](#)

Ordering Information

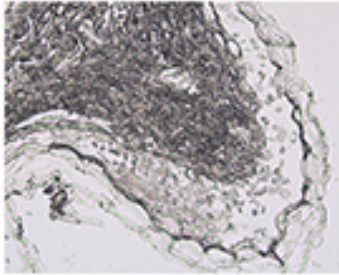
[Order Online »](#)

ENZ-C34904	~10 µg - 0.5ml
------------	----------------

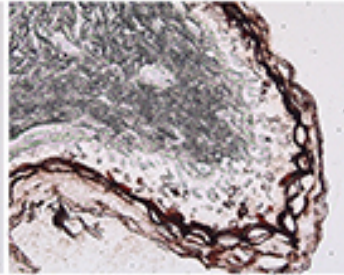
Manuals, SDS & CofA

[View Online »](#)

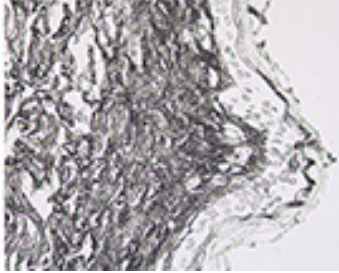
(-) control, 1X Dilution Buffer



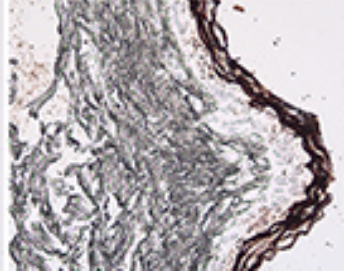
(+) ENZ-C34904



(-) control, 1X Dilution Buffer



(+) ENZ-C34904



Immunohistochemistry analysis of human skin tissue using Cytokeratin 1 (human), mAb (34 β B4) at a dilution of 1:20.

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CK1
Application	IHC (PS)
Clone	34βB4
Formulation	Liquid. 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). A 20- to 50-fold dilution of the antibody is suggested. 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 .



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)