

Cytokeratin 14 polyclonal antibody

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

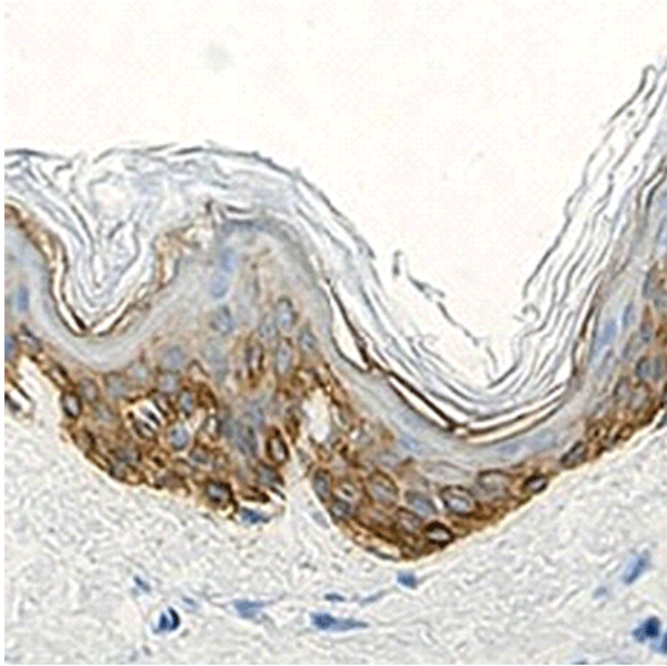
Ordering Information

[Order Online »](#)

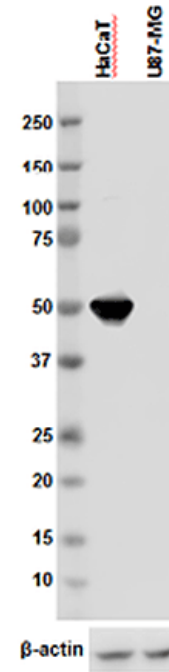
ENZ-ABS171A-0100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



IHC staining of anti-Keratin 14 antibody (Poly19053) on formalin-fixed paraffin embedded normal human skin.



Western blot analysis of cell lysates from HaCaT (Positive control) and U87-MG (Low expression negative control) cells using Keratin 14 Rabbit primary antibody and HRP Donkey anti-Rabbit secondary antibody.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name CK14

Application ICC, IHC, WB

Crossreactivity Dog, Primate

Formulation Liquid. In phosphate-buffered solution, pH 7.2, containing 0.09% sodium azide.

Host Rabbit

Immunogen Peptide sequence derived from the C-terminus of mouse keratin 14.

Isotype IgG

Purity Detail Affinity purified.

Species Reactivity Human, Mouse, Rat

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)