

Filaggrin polyclonal antibody (ENZ-ABS181)

IF

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Skin	Keratinocyte cell line	HaCaT	https://www.ncbi.nlm.nih.gov/pubmed/29247241
Mouse	Skin	Dorsal skin tissue section		https://www.ncbi.nlm.nih.gov/pubmed/30132547

For Technical Support:
techserv-usa@enzolifesciences.com
techserv-eu@enzolifesciences.com

Complete citation can be found on product's webpage

Global Headquarters
ENZO LIFE SCIENCES, INC.
Ph: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES (ELS) AG
Ph: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Ph: +32 3 466 0420
info-be@enzolifesciences.com

France
Ph: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Ph: +49 7621 5500 526
info-de@enzolifesciences.com

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

For local distributors and detailed product information visit us online: www.enzolifesciences.com

For Research Only. Not for Use in Diagnostic Procedures.