

Gelsolin (human) monoclonal antibody (20)

Gelsolin participates in the extracellular actin scavenger system, binding actin and disaggregating filamentous actin released from dying cells. Actin monomers are then transferred to Gc-globulin. The plasma concentration of gelsolin is reduced in conditions of major actin release, such as acute liver failure, major trauma and rhabdomyolysis, where multi-organ failure is a risk. It has been suggested that gelsolin, like Gc-globulin, can be a prognostic factor in these conditions.

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

Manuals, SDS & CofA

[View Online »](#)

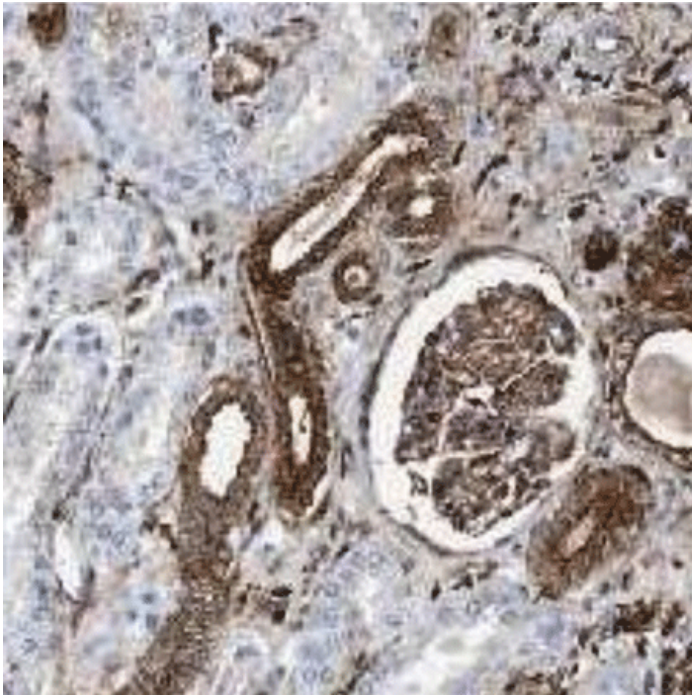


Figure 2. Immunohistochemical staining of human kidney with ABS 017-20 shows cytoplasmic and membranous positivity in both renal tubules and glomeruli. Antibody dilution: 1:6000

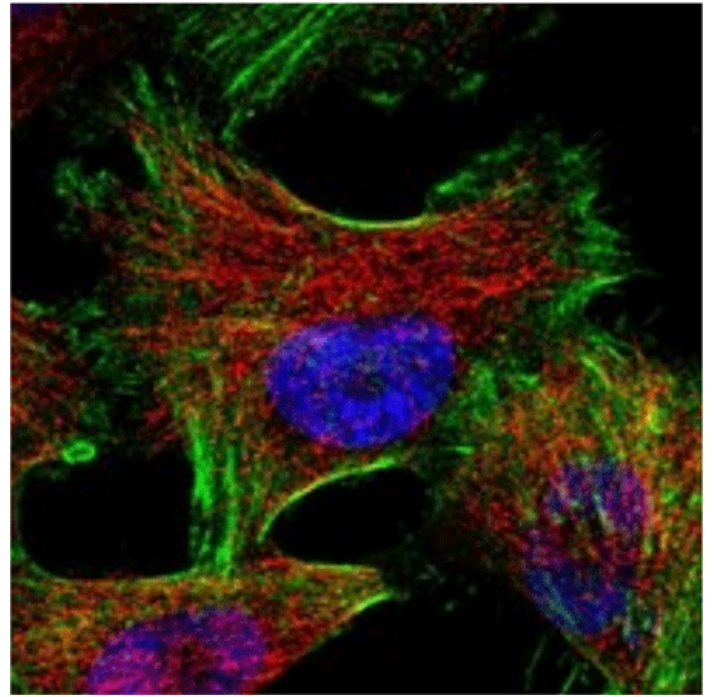


Figure 1. Immunofluorescent staining of human cell line U-251 MG with BPD-ABS-017-20 shows positivity in actin filaments. Antibody dilution: 1:2000

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	ELISA, IF, IHC, WB
Clone	20
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.
GenBank ID	2934
Host	Mouse
Immunogen	Purified gelsolin from human plasma.
Isotype	IgG1κ
Purity Detail	Protein A/G-purified.
Source	From culture supernatant.
Species Reactivity	Human
UniProt ID	P06396
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

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

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

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

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

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