

MET monoclonal antibody (BFA-13)

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

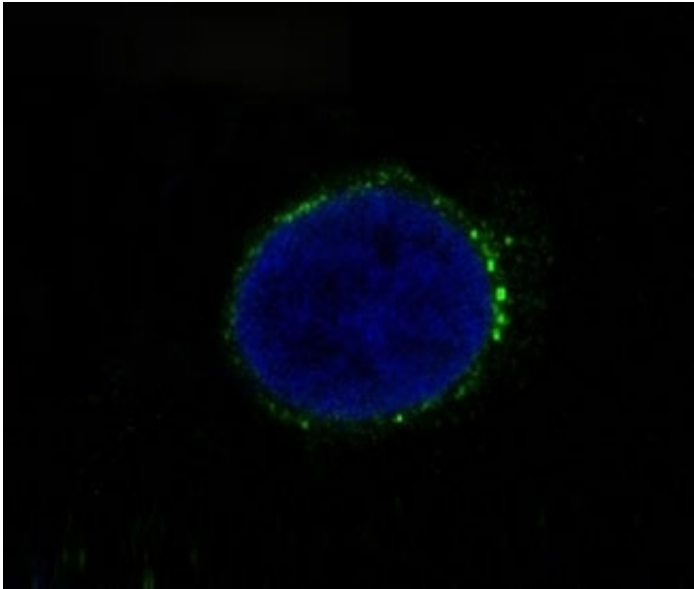
Ordering Information

[Order Online »](#)

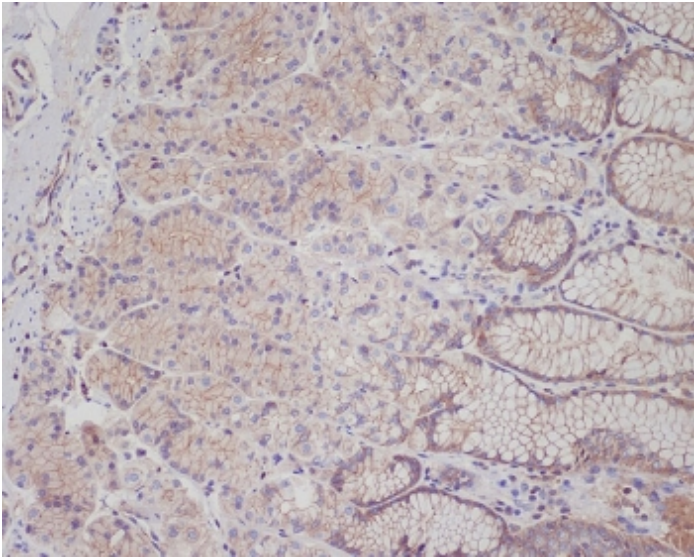
ENZ-ABS905-0100	100µl
-----------------	-------

Manuals, SDS & CofA

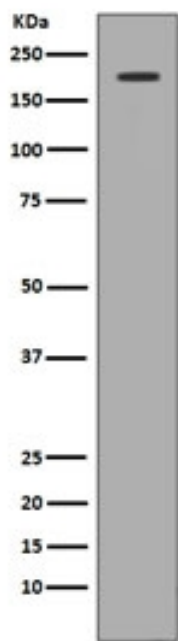
[View Online »](#)



Immunofluorescent staining of FFPE human HT-29 cells with c-Met antibody (green) and DAPI nuclear stain (blue). HIER: steam section in pH6 citrate buffer for 20 min.



IHC staining of FFPE human stomach tissue with c-Met antibody. HIER: boil tissue sections in pH6, 10mM citrate buffer, for 10-20 min and allow to cool before testing.



Western blot testing of human 293 cell lysate with c-Met antibody. Predicted molecular weight ~156 kDa.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	c-Met, HGFR, Hepatocyte growth factor receptor
Application	IF, IHC (PS), WB
Clone	BFA-13
Formulation	Liquid. In PBS containing 0.02% sodium azide., 50% glycerol and 0.4-0.5mg/ml BSA
Host	Rabbit
Immunogen	Synthetic peptide specific to human c-Met / MET.
Isotype	IgG
Purity Detail	Affinity purified.
Recommendation Dilutions/Conditions	Immunofluorescence (1:50-1:200)Immunohistochemistry (FFPE) (1:50-1:200)Western blot (1:500-1:2,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	P08581
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee



Last modified: July 12, 2024

ENZO LIFE SCIENCES,
INC.
Phone: +1 940 945 1139
info@enzolifesciences.com
usa@enzolifesciences.com

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

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

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

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

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