

Birch Pollen monoclonal antibody (202B8.03) (purified)

Ordering Information

[Order Online »](#)

DEN-DDX0052P-100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Regulatory Status RUO - Research Use Only

Product Details

Application Notes	Binding to nitrocellulose-blotted Bet v 1, inhibition of human IgE binding to Bet v1, cross-competition with anti-Bet v1 mAbs, Bet v1 immunoprecipitation
Clone	202B8.03
Host	Mouse
Isotype	IgG1
Species Reactivity	Birch
Technical Info / Product Notes	For the Original Manufacturer's data sheet please click here .



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