

Bordetella pertussis

FHA

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-630-123-0100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

- Minimal degradation following extraction

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Filamentous hemagglutinin antigen
Appearance	White powder.
Formulation	Lyophilized from 0.05M TRIS-HCl and 0.5M sodium chloride at pH 8, 5% sucrose.
Purity	≥98% (SDS-PAGE)
Source	From <i>Bordetella pertussis</i> , strain NCTC 10739.
UniProt ID	P12255

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