

Cryo-EM Puck Basic Starter Kit

Ordering Information

[Order Online »](#)

JBS-X-CEMG2-BLSK	1Kit
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Contents

- 10 Cryo-EM Pucks Gen 2.0 (X-CEMG2-10)
- 1 Cryo-EM Puck Storage Cane for HC34 & VHC35 (X-CEM-202)
- 1 Cryo-EM Puck Shipping Cane for CX100 (X-CEM-204)
- 1 Cryo-EM Puck Grasping Tongs (X-CEM-GT)



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