

Cryo-EM Puck Gen 2.0

(10 pcs.)

The Cryo-EM Pucks Gen 2.0 have enhanced features over other cryo-EM pucks on the market. These include: 12 numbered well locations with easy-access tweezer notches Magnets in the bottom of the pucks, which securely hold the puck to the shelves of the shipping and storage canes Front grabbing tong location 2D barcoding for advanced sample tracking

Ordering Information

[Order Online »](#)

| | |
|----------------|------|
| JBS-X-CEMG2-10 | 10ea |
|----------------|------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details



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