

XtalTool HT 18 mm

Sample Holder for Crystal Growth, in situ Ligand Soaking and in situ Data Collection with a Robot Assisted Sample Mounting

The XtalTool HT is a patent-pending sample holder consisting of a plastic support with an outer transparent COC film and an inner X-ray transparent yellow polyimide film with 5 μm pores. The design allows the setup of two crystallization drops on one XtalTool HT. Crystal growth can be monitored either by a transmission microscope or an automated inspection system. All intended manipulative steps on the crystal are carried out in-situ. Before data collection the outer ring can easily be broken off at the designated respective break points. Mounted on an XtalTool Base, the XtalTool HT then fits into a SPINE standard magnetic CryoVial and thereby facilitates robot assisted sample mounting. Its geometry meets the 18 mm SPINE standard and it's compatible with most synchrotron and in-house beamlines dedicated to macromolecular crystallography.

Ordering Information

[Order Online »](#)

JBS-X-XT-104	20ea
--------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

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