# **XtalTool** The ALL-in-ONE Sample Holder





Crystal handling is a major bottleneck in protein crystallography. Macromolecular crystals are often fragile and suffer from mechanical stress during manipulation and/or mounting. This can lead to impairment and even complete loss of diffraction quality. **XtalTool** was developed to minimize the handling steps of crystals in order to obtain the best possible diffraction data<sup>[1]</sup>.

### 1. XtalTool is used as a cover slide for hanging drop crystallization









2. Crystal manipulation such as ligand soaking or cryoprotection is performed in a gentle way directly on the sample holder



#### 3. X-ray diffraction data are collected in-situ at ambient or cryo temperature



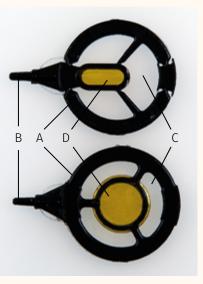
# **XtalTool** The ALL-in-ONE Sample Holder





## One technology - two designs

The patent-pending XtalTool consists of a round black plastic support (A) with an outer diameter of 18 or 22 mm and an extension (B) that fits into a customized goniometer base. A transparent removable COC film (C) tightly seals the well during crystallization. An inner X-ray transparent polyimide film (yellow) with 5 µm pores (D) enables a gentle crystal soaking procedure by diffusion and serves as crystal support during data collection.



**XtalTool HT** with predetermined breaking points at the outer frame is compatible with robot assisted sample mounting. It is available in 22 mm diameter (Cat.-No. X-XT-103) to fit the Crystalgen SuperClear Plates and in 18 mm diameter (Cat.-No. X-XT-104) to fit the smaller Greiner Combo Plates with SBS footprint.

The classic **XtalTool** with 22 mm diameter (Cat.-No. X-XT-101) can be used with Crystalgen SuperClear Plates and is manually mounted on the goniometer.

Product	CatNo.	Amount
XtalTool HT 22 mm	X-XT-103	20 pcs.
XtalTool HT 18 mm	X-XT-104	20 pcs.
XtalTool Bases	X-XT-105	20 pcs.
XtalTool	X-XT-101	24 pcs. + 1 Base
XtalTool Soaking Kit	X-XT-102	1 Kit

XtalTool is unique on the market, it was developed by the HZB MX-group at BESSY II (AG Weiss).

#### **Reference:**

[1] Feiler et al. (2019) An All-in-one Sample Holder for Macromolecular X-ray Crystallography with Minimal Background Scattering. J. Vis. Exp. 149:e59722.



Jena Bioscience GmbH Loebstedter Str. 71 07749 Jena Germany Phone +49(0)3641-6285 000 Fax +49(0)3641-6285 100 info@jenabioscience.com www.jenabioscience.com



Certified QMS and EMS according to DIN EN ISO 9001 and DIN EN ISO 14001 Reg.-No.: ICV03597 034 and ICV03597 534

Crystallography & Cryo-EM