

**Quantifoil™ Multi Holes on 300 copper mesh**

Cat. No.	Amount
X-160-CU300	100 pcs.

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a

Applications:

Support film for cryo-tomography (cryo-ET) or cryo-electron microscopy (cryo-EM).

Description:

QUANTIFOIL™ Multi Holes offer a pattern of various hole sizes, shapes and arrangements applied on 300 copper mesh. The regular repetition allows the use of modern automated data collection strategies.

Hole dimensions:

- Circular holes have diameters of 1 μm , 1.4 μm , and 2 μm
- Elliptical holes measure 1 μm x 4 μm and 2 μm x 8 μm , they appear round in tomographic sections with high tilt angles of 70°