

**Quantifoil™ R 1.2/1.3 on 300 copper mesh**

Cat. No.	Amount
X-101-Cu300	100 pcs.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a

Applications:

Support film for biological samples in Cryo-EM techniques

Description:

QUANTIFOIL™ R 1.2/1.3 carbon film has circular holes of 1.2 µm diameter with an interspace of 1.3 µm in both dimensions, applied on 300 copper mesh. Suitable for magnifications around 50.000x and higher.