

**Quantifoil™ R 2/2 on 300 copper mesh**

Cat. No.	Amount
X-103-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 biological samples in Cryo-EM techniques

Description:

QUANTIFOIL™ R 2/2 carbon film has circular holes of 2 µm diameter with an interspace of 2 µm in both dimensions, applied on 300 copper mesh. Suitable for magnifications between 30.000x and 40.000x.