

**Quantifoil™ R 3.5/1 on 200 copper mesh**

Cat. No.	Amount
X-104-Cu200	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 3.5/1 carbon film has circular holes of 3.5 µm diameter with an interspace of 1 µm in both dimensions, applied on 200 copper mesh. The larger hole size makes these grids useful for imaging cells and viruses, they are suitable for magnifications around 20.000x.