

**Quantifoil™ R 3.5/1 on 300 copper mesh**

Cat. No.	Amount
X-104-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 3.5/1 carbon film has circular holes of 3.5 µm diameter with an interspace of 1 µm in both dimensions, applied on 300 copper mesh. The larger hole size makes these grids useful for imaging cells and viruses, they are suitable for magnifications around 20.000x.