

**Quantifoil™ R 1/4 on 200 copper mesh**

Cat. No.	Amount
X-105-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 1/4 carbon film has circular holes of 1 µm diameter with an interspace of 4 µm in both dimensions, applied on 200 copper mesh.