

**Quantifoil™ R 1.2/1.3 plus C2 on 300 copper mesh**

Cat. No.	Amount
X-101-Cu300C2	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 plus C2 is a 12 nm carbon film with circular holes of 1.2 µm diameter and an interspace of 1.3 µm in both dimensions, which has an additional 2 nm continuous carbon layer on top. Applied on 300 copper mesh.