

**Quantifoil™ R 2/1 plus C2 on 300 copper mesh**

Cat. No.	Amount
X-102-Cu300C2	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 2/1 plus C2 is a 12 nm carbon film with circular holes of 2 µm diameter and an interspace of 1 µm in both dimensions, which has an additional 2 nm continuous carbon layer on top. Applied on 300 copper mesh.