## **DATA SHEET**





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

Cat. No.	Amount
X-101-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<sup>TM</sup> R 1.2/1.3 plus C2 is a 12 nm carbon film with circular holes of 1.2  $\mu$ m diameter and an interspace of 1.3  $\mu$ m in both dimensions, which has an additional 2 nm continuous carbon layer on top. Applied on 300 copper mesh.