

**Quantifoil™ R 2/1 on 300 gold mesh**

| Cat. No.    | Amount   |
|-------------|----------|
| X-102-Au300 | 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 2/1 carbon film has circular holes of 2 µm diameter with an interspace of 1 µm in both dimensions, applied on biocompatible 300 gold mesh. Suitable for magnifications between 30.000x and 40.000x.