



## C-flat™ 2/1 on 300 gold mesh

Cat. No.	Amount
X-302-AU300	50 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:**

20 nm C-flat™ carbon film with 2 µm hole size and 1 µm hole spacing, applied on 300 gold mesh.

**nonsaleable to USA**