



C-flat™ 2/1 on 400 gold mesh

Cat. No.	Amount
X-302-AU400	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 400 gold mesh.

nonsaleable to USA