

Proteomics

Using the proteoCHIP 12*16 autosampler holder for the Vanquish Neo UHPLC system

Authors

Runsheng Zheng¹, Bernard Delanghe²,
Christopher Pynn¹, David Hartlmayr³,
Anjali Seth³

¹Thermo Fisher Scientific,
Germering, Germany

²Thermo Fisher Scientific,
Bremen, Germany

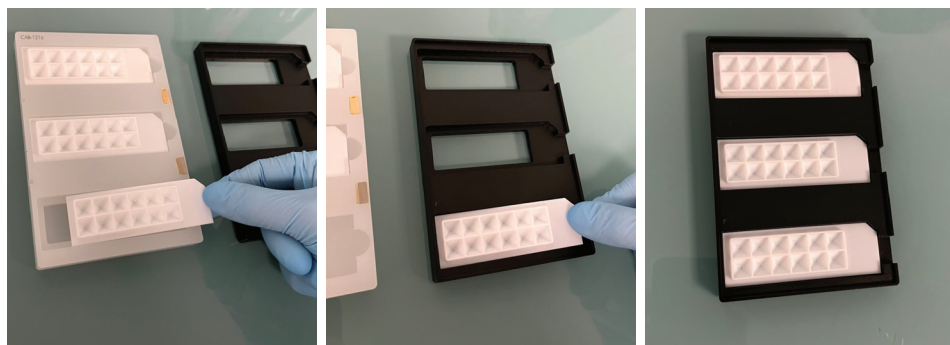
³Cellenion, Lyon, France

Keywords

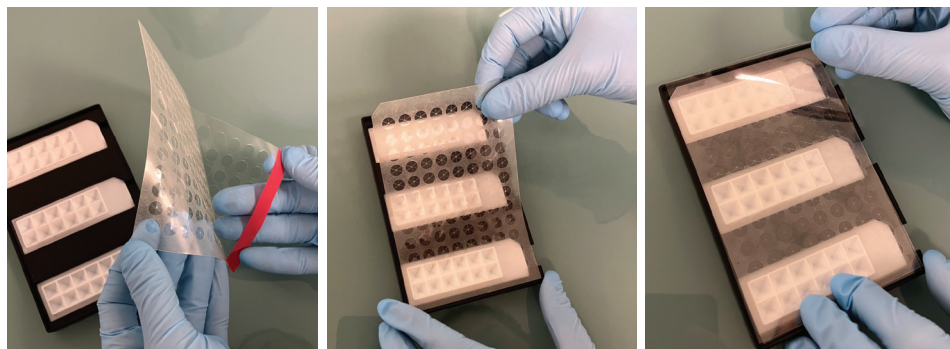
Orbitrap mass spectrometer,
Vanquish Neo UHPLC system,
single-cell proteomics, cellenONE,
TMTpro 16plex, proteoCHIP, bottom
detection, SCoPE-MS

The Cellenion proteoCHIP 12*16 autosampler holder is used to interface the proteoCHIP 12*16 funnels with the Thermo Scientific™ Vanquish™ Neo UHPLC system for Thermo Scientific™ TMTpro™ 16plex label-based LC-MS analysis. The following procedure details its use.

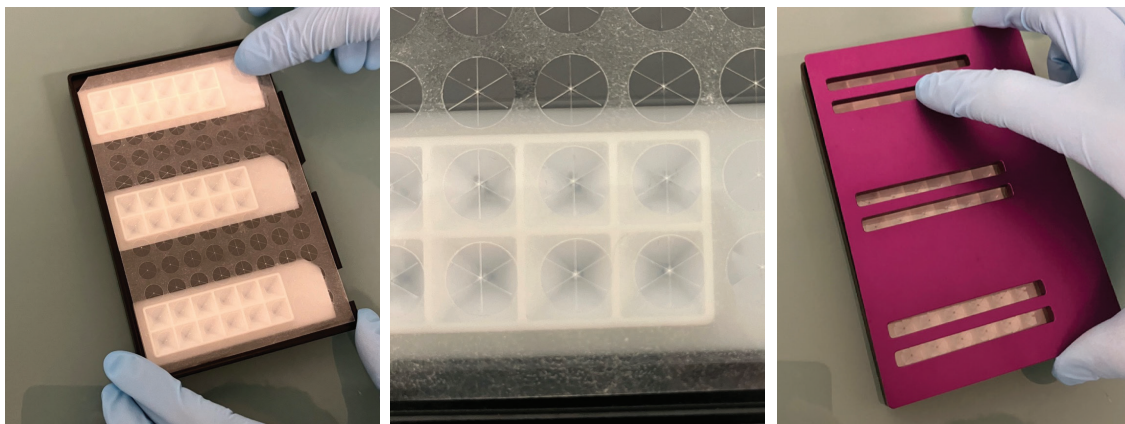
1. After centrifugation of the TMTpro-labeled samples from the proteoCHIP 12*16, transfer proteoCHIP 12*16 funnels containing the pooled samples to the proteoCHIP 12*16 autosampler holder base (black).



2. Remove the bottom protective film (on the side with the adhesive) of the sealing tape (Thermo Fisher Scientific, P/N 60180-M146/M176). Position the sealing tape over the funnels and peel off the upper protective film.

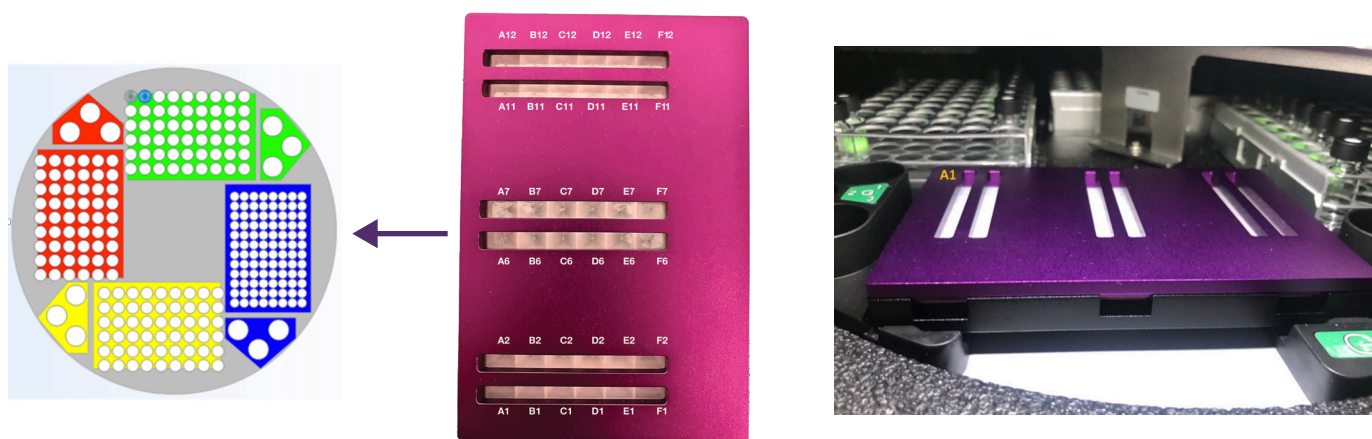


3. Make sure that the pre-slit circles of the sealing tape are well aligned to the center of the funnels and place the holder lid (pink) on top.



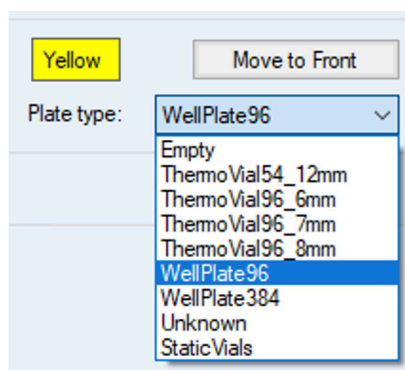
4. Position the holder in the autosampler carousel of the Vanquish Neo UHPLC system such that the “A1” label is in the top left-hand corner.

Warning: Failure to insert the adapter in the correct orientation could result in damage to the autosampler needle.



5. To enable the sampler to recognize the holder for sample injection, set the plate type to “WellPlate96” manually on the SII e-panel in the MS controlling PC.

Warning: Ensure that this setting is applied to all sampler trays where the proteoCHIP 12*16 autosampler holder is present.



6. Inject samples **exclusively from the specified positions listed below** during LC-MS analysis. There are no additional sampler settings required on the LC.

Left funnel: A1, B1, C1, D1, E1, F1; A2, B2, C2, D2, E2, F2

Middle funnel: A6, B6, C6, D6, E6, F6; A7, B7, C7, D7, E7, F7

Right funnel: A11, B11, C11, D11, E11, F11; A12, B12, C12, D12, E12, F12



Warning: Failure to adhere to these positions may result in damage to the LC sampler needle.

Learn more at thermofisher.com/singlecellproteomics