

August 19, 2020

To Whom it may concern,

In early 2019, we implemented a new pump configuration for the 42i instruments for improved quality and performance.

The new pump retrofit kit, PN 119775-00, replaces the legacy internal pump, PN 101011-00 and 101011-01 in all 42i configurations. The external pump in some options remains unchanged.

Improvements include:

- Stainless steel valve plate – more consistent quality over Teflon valve plate
- PTFE Impregnated Hard Anodized Intermediate plate – better sealing for deeper vacuum

If you have any questions, please feel free to contact me.

Sincerely,



Joseph Vaglio

Product Manager, Gas Analyzers

Mobile: +1 (508) 954-1715

Joseph.Vaglio@thermofisher.com