



June 9, 2015
UVPM: 43,098

**Quanterix Multiplex Simoa Assays and Homebrew Kit
By Andrew Han**

Quanterix announced the launch of its Simoa-based multiplex panels and homebrew kit. The first commercially available multiplex kits are 3-plex panels of cytokines for measuring inflammatory response across a wide range of conditions, while the homebrew kit allows customers to use their own antibody pairs to develop custom multiplex Simoa panels.