

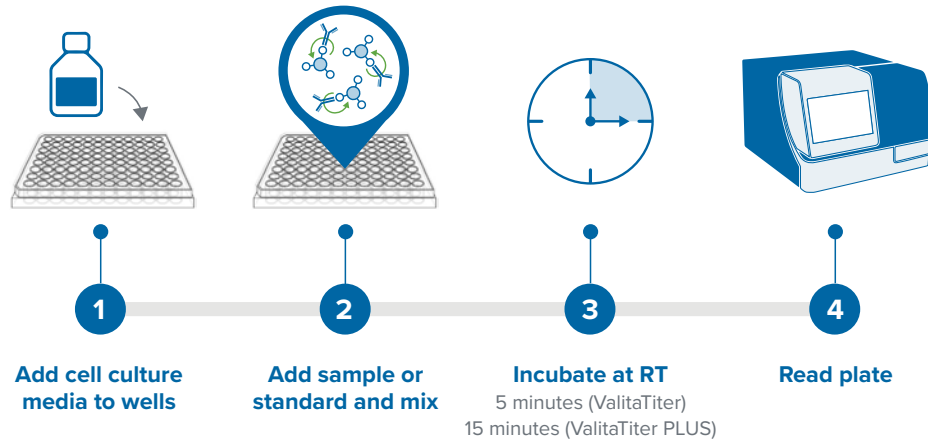
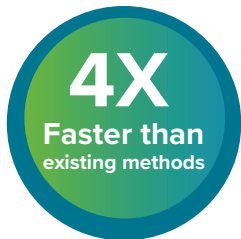


Quickly and easily quantify IgG with ValitaTiter

Integrating ValitaTiter and Molecular Devices fluorescence polarization (FP) plate readers produces a seamless IgG titer assay setup and optimization. Cell line development teams will realize savings on time and resources, as well as faster time-to-market by utilizing a single technology for measuring a critical quality attribute during target selection, process development, and production QC release in (automated) high-throughput workflows.

ValitaTiter workflow

Standard curve workflow:



Support

Molecular Devices' technical support team and field application scientists elevate the ValitaTiter customer experience with deep knowledge of protocol optimizations for ValitaTiter applications.



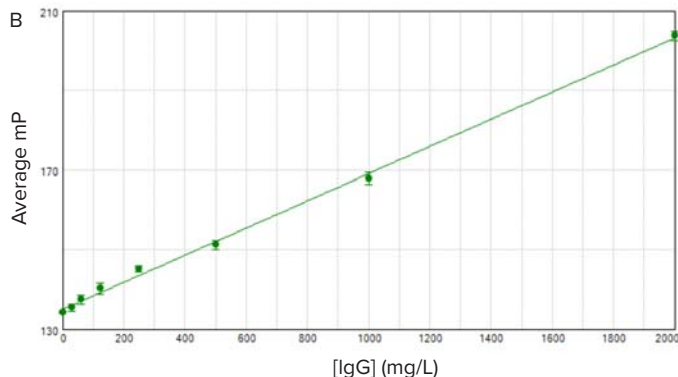
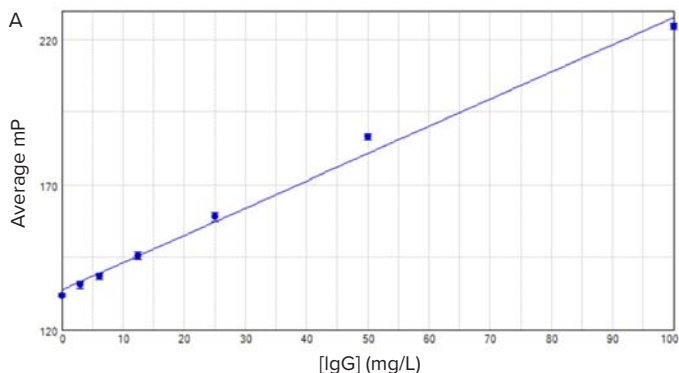
Automation

In-house automation experts provide ValitaTiter customers robotic workflow customization and automation.



GxP

GxP compliance software ensures regulatory guidelines are met in a manufacturing site.



Standard curves for the ValitaTiter (A, $r^2 = 0.993$) and ValitaTiter Plus (B, $r^2 = 0.998$) assays. Curves were plotted using a linear curve fit in SoftMax[®] Pro Software. *Quantifiable IgG ranges for kit: ValitaTiter: 2.5 mg/L -100 mg/L; ValitaTiter Plus: 100 mg/L - 2000 mg/L

Compatible microplate readers

	iD5	i3x	M5	Paradigm
Delta mP	93	70	80	89
Average StDev (mP)	1.12	2.12	0.57	0.82
Average %CV	0.7	0.7	0.5	0.3
Max StDev	1.54	3.53	1.06	1.27
Max %CV	1.0	1.3	1.0	0.4

Standard delta mP, average standard deviation (StDev) and %CV (n = 4) for ValitaTiter standards read on Molecular Devices readers.

	iD5	i3x	M5	Paradigm
Delta mP	69	85	92	88
Average StDev (mP)	1.05	3.54	1.63	1.14
Average %CV	0.7	1.3	1.6	0.4
Max StDev	1.59	6.49	2.26	1.43
Max %CV	1.1	2.0	2.4	0.5

Standard delta mP, average standard deviation (StDev) and %CV (n = 4) for ValitaTiter Plus standards read on Molecular Devices readers.

See our app note for details: <https://www.moleculardevices.com/en/assets/app-note/br/high-throughput-igg-quantitation-platform-for-clone-screening-during-drug-discovery-and-development>

ValitaTiter comparison

	ValitaTiter	BLI	ELISA	HPLC
Assay time	15 minutes	1 hour	5 hours	10 hours
Sample prep	None	None	Cell centrifugation	Cell centrifugation and protein purification
Reagents	1	>2	>6	2
Steps	Add, mix, read	>3	>20 steps	>5
Costs	\$	\$\$\$\$	\$	\$\$\$

Ordering information

Product Name	Part Number
ValitaTiter 96-well (1 plate)	VAL003
ValitaTiter Plus 96-well (1 plate)	VAL004
ValitaTiter 96-well (Pack of 20)	VAL003 PACK
ValitaTiter Plus 96-well (Pack of 20)	VAL004 PACK

Phone: +1.877.589.2214
 Web: www.moleculardevices.com
 Email: NSD@moldev.com
 Check our website for a current listing of worldwide distributors.

USA and Canada +1.800.635.5577
 United Kingdom +44.118.944.8000
 Europe* 00800.665.32860
 China +86.4008203586

*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Luxembourg, Netherlands, Portugal, Spain, Sweden, Switzerland and United Kingdom

Compatible Fluorescence Polarization Plate Readers

FP Plate Reader
SpectraMax [®] M5/M5e Multi-Mode Microplate Reader
SpectraMax [®] iD5 Multi-Mode Microplate Reader
SpectraMax [®] i3x Multi-Mode Microplate Reader
SpectraMax [®] Paradigm Multi-Mode Microplate Reader
SpectraMax [®] FlexStation 3 Multi-Mode Microplate Reader

Taiwan/Hong Kong +886.2.2656.7585
 Japan +81.3.6362.9109
 South Korea +82.2.3471.9531
 India +91.73.8661.1198