

Capability/Features	Monolith X	Monolith Omni
In-solution affinity measurement	✓	✓
Spectral Shift	✓	✓
TRIC	✓	✓
Kinetics rates (k _{on} , k _{off})	—	✓
Ligand-induced stability analysis	—	✓
Thermodynamic parameters (ΔH, ΔG)	✓	✓
Best fit	Low-sample affinity measurements when K _d is the primary decision point	Complete interaction profiling beyond K _d , with affinity, kinetic rates, and ligand-induced stability insight

Talk to a specialist

Find the right configuration for your needs

Contact Us →