



SpectraMax[®] i3x

Multi-Mode Microplate Reader

A flexible and reliable high-performance microplate reader for your virus research.

Nucleic acid quantification

Measure nucleic acid concentration directly using absorbance; or use fluorescence for greater sensitivity

Cell signaling assays

Assess cellular response to viral infection

Biomolecular binding

Study viral assembly, other viral processes; detect viral components, viral inhibition

Total protein quantification (non-specific)

Measure protein (e.g., viral protein) concentration directly using absorbance; or use a colorimetric method for greater sensitivity

Assays to quantify viruses and study viral processes

Measure viral replication, virus neutralization

Cell viability

Assess cell viability by measuring ATP levels or cellular enzyme activity

Detection and quantification of specific proteins

Detect and measure viral antigens, or immunoglobulin produced in response to virus exposure

[Click here to contact your local sales representative to find out more.](#)

Committed to supporting scientists in their research of COVID-19 cellular response and vaccine development.

Contact Us

Phone: +1.800.635.5577
 Web: www.moleculardevices.com
 Email: info@moldev.com

Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada	+1.800.635.5577	China (Beijing)	+86.10.6410.8669	Japan	+81.3.6362.9109
United Kingdom	+44.118.944.8000	China (Shanghai)	+86.21.3372.1088	South Korea	+82.2.3471.9531
Europe*	00800.665.32860	Hong Kong	+852.3971.3530		

*Austria, Belgium, Denmark, Finland, France, Germany, Ireland, Netherlands, Spain, Sweden and Switzerland