



revvity

Detect mpox with confidence.

Mpox Virus Real Time PCR Kit

Our kit is designed to detect the mpox virus, including the new strain responsible for the recent outbreak. With highly selective targeting, it focuses on a stable region of the mpox genome, ensuring reliable virus detection.

Our kit provides:



Reliable detection: Our kit includes both positive and negative internal controls for reliable analysis of results and detects both clade I and II mpox virus, increasing confidence in the result



Highly selective targeting The test targets a highly conserved region of the mpox genome (F3L), resulting in reliable analysis



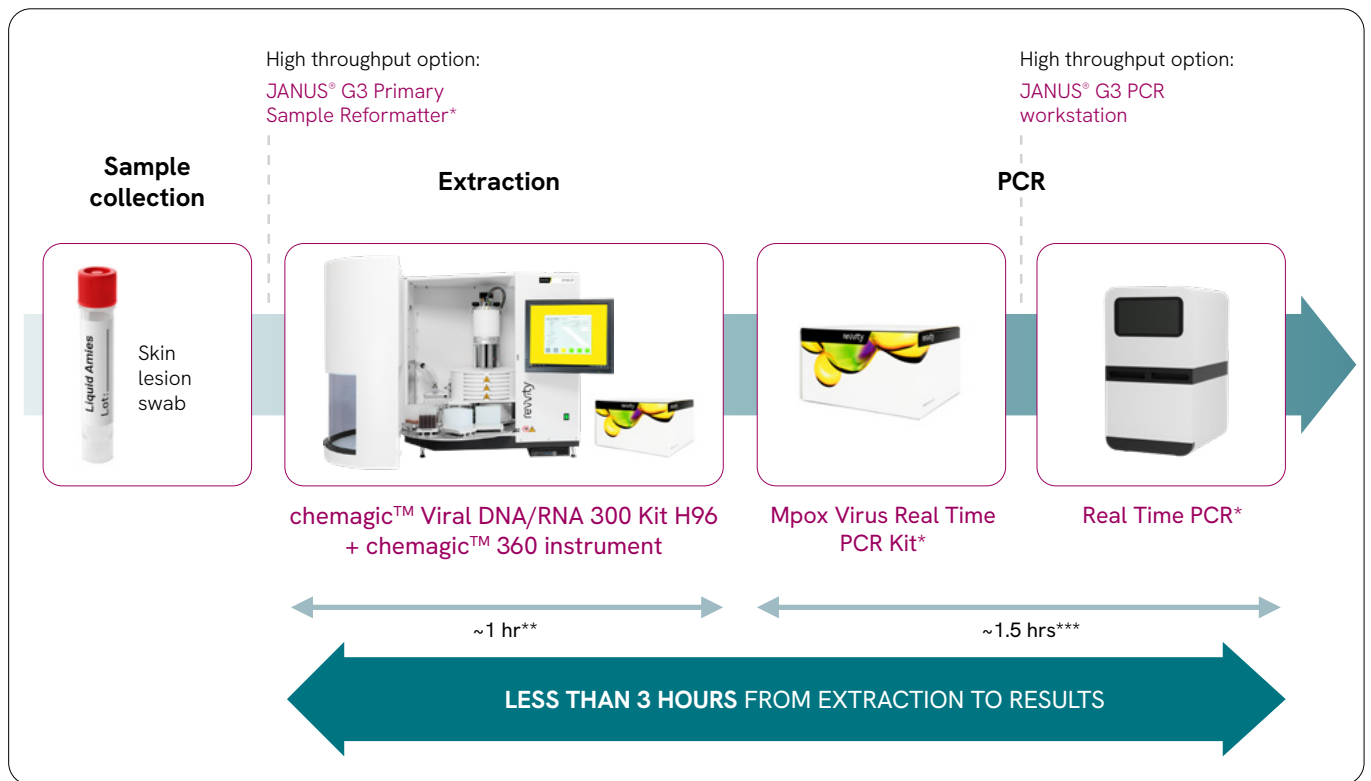
Fast results: With a quick turnaround time of less than 3 hours from sample extraction to results, our workflow effectively supports essential research in a timely manner

Contact us today:



Automated workflows

With Revvity's automation solutions, you can receive your mpox results in **less than 3 hours** from sample extraction to analysis.



* This assay has been validated on commonly used PCR instruments.

** Validated extraction protocol of 1 hour on a full 96 well plate.

*** Based on full run using one plate of 96 samples.



Extensive expertise and capacity

With our extensive expertise in infectious diseases and our large manufacturing capabilities, we can quickly develop and deliver products efficiently to address global emergencies.

revvity