

Luminex



xTAG Respiratory Viral Panel FAST

Actionable results for better patient care*

96 samples in 5 hours

The xTAG[®] RVP FAST is a qualitative molecular diagnostic test for the simultaneous detection of multiple viral types and subtypes from nasopharyngeal swabs of individuals suspected of respiratory tract infections.

Streamlined Workflow

- Reduced hands on time and faster workflow allow you to perform comprehensive front-line respiratory testing
- Consolidated testing of multiple respiratory viruses into a single assay improves labor efficiency and cost-effectiveness
- Scalable assay built on the versatile Luminex[®] 100/200™ platform

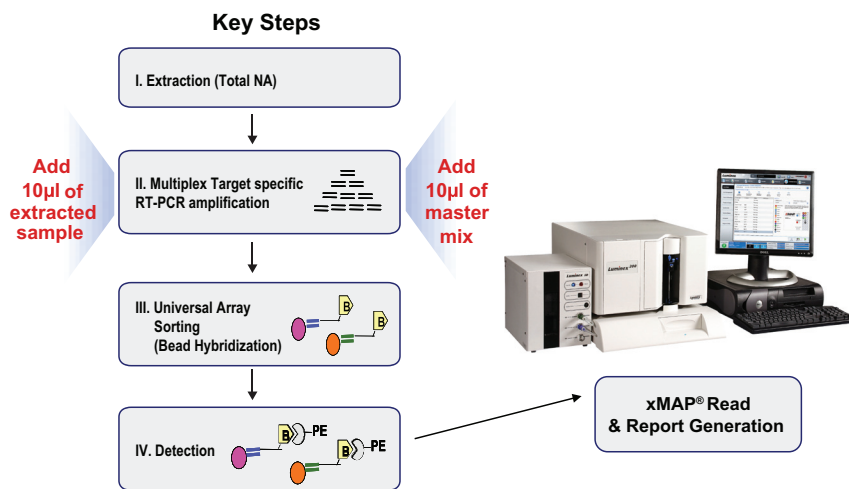
Actionable Results*

- More comprehensive than most PCR based front-line tests
- Confidence in performance with clinical sensitivity and specificity that exceeds 90% for viral targets

Viral Types and Subtypes detected by RVP FAST

Respiratory Syncytial Virus (RSV)
Influenza A
Non-specific influenza A
H1 subtype
H3 subtype
Influenza B
Human Metapneumovirus (hMPV)
Rhinovirus
Adenovirus
<i>MS-2 Bacteriophage Internal Control</i>
<i>Bacteriophage Lambda Positive Control</i>

Assay Workflow



The xTAG RVP FAST Assay is distributed in the US from:

Abbott Molecular Inc. Catalog No. 5N01-01 Kit size: 96 tests	Fisher HealthCare Catalog No. 22-575-470 Kit size: 96 tests	Luminex Catalog No. I025C0197 Kit size: 96 tests
--	---	--

For More Information:

Visit the Luminex website at www.luminexcorp.com
Call 1-800-593-2370 or email info@luminexcorp.com
for sales and product information.

IVD For In Vitro Diagnostic Use

UNITED STATES
Luminex Corporation
Austin, Texas
Tel: 512.219.8020
Fax: 512.219.5195

CANADA
Luminex Molecular Diagnostics
Toronto, Ontario
Tel: 416.593.4323
Fax: 416.593.1066

EUROPE
Luminex B.V.
Oosterhout, The Netherlands
Tel: +31.162.408333
Fax: +31.162.408337

CHINA
Luminex Shanghai Trading Co.
Shanghai, China
Tel: +86.21.616.50809
Fax: +86.21.616.50811

JAPAN
Luminex Japan Corporation, Ltd.
Tokyo, Japan
Tel: +81.3.5545.7440
Fax: +81.3.5545.0451

AUSTRALIA
BSD Robotics, A Luminex Company
Brisbane, Queensland
Tel: +61.7.3273.0273
Fax: +61.7.3273.0274

www.luminexcorp.com

info@luminexcorp.com