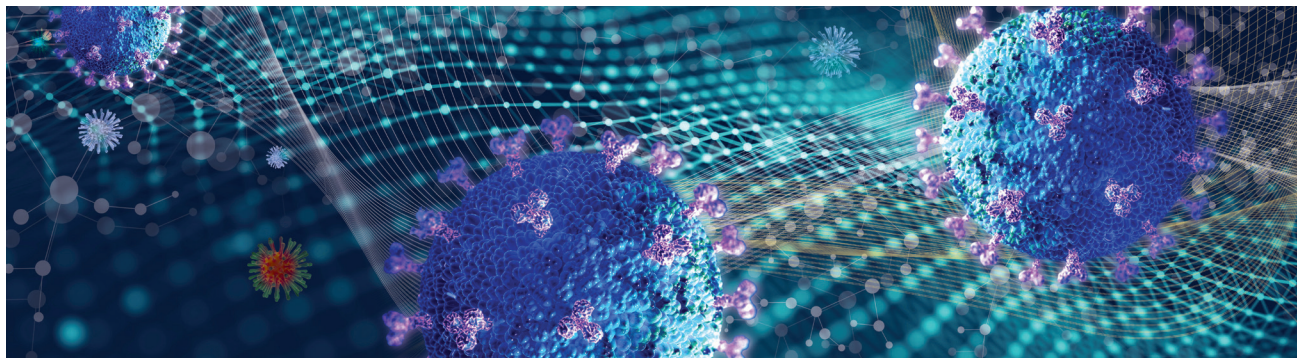




Bio-Rad Launches the SARS CoV-2 Standard for Coronavirus (COVID-19) Research Testing

The SARS CoV-2 Standard and SARS CoV-2 Negative from the Exact Diagnostics product line contain genomic DNA allowing laboratories to validate the entire process of a molecular assay including extraction, amplification, and detection.

SARS-CoV-2 Standard

The EDX SARS-CoV-2 Standard is manufactured with synthetic RNA transcripts containing 5 gene targets (E, N, S, ORF1a, and RdRP genes of SARS-CoV-2, 200,000 copies/mL each). The product is formulated in a synthetic matrix and contains human genomic DNA (75,000 copies/mL).

SKU: COV019

5 Vials, 0.3 mL Fill Volume, -20°C or below Storage



SARS-CoV-2 Negative

The EDX SARS-CoV-2 Negative is negative for SARS-CoV-2. The product is formulated in a synthetic matrix and contains human genomic DNA (75,000 copies/mL)

SKU: COV000

5 Vials, 0.3 mL Fill Volume, -20°C or below Storage



LIMITATIONS: For research use only. Not for use in diagnostic procedures.



Contact your local Bio-Rad sales representative for more information