





SKU/PRODUCT NAME	DESCRIPTION	INTENDED USE
TRANSPLANTATION		
ADVP100 Adenovirus Verification Panel	5 Panel Members, 1.0 mL, -20°C Storage, Dynamic Range 200 cp/mL - 2,000,000 cp/mL	CE-IVD, IVD
ADVL101 Adenovirus Low Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	CE-IVD, IVD
ADVH102 Adenovirus High Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	CE-IVD, IVD
BKVP100 BKV Verification Panel	6 Panel Members, 1.0 mL, -20°C Storage, Dynamic Range 200 IU/mL - 20,000,000 IU/mL	CE-IVD, IVD
BKVL101 BKV Low Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	CE-IVD, IVD
BKVH102 BKV High Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	CE-IVD, IVD
B19P100 Parvovirus B19 Verification Panel	5 Panel Members, 1.0 mL, -20°C Storage, Dynamic Range 500 IU/mL - 5,000,000 IU/mL	CE-IVD, IVD
B19L101 Parvovirus B19 Low Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	CE-IVD, IVD
B19H102 Parvovirus B19 High Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	CE-IVD, IVD
CMVP100 CMV Verification Panel 1.0 mL	5 Panel Members, 1.0 mL, -20°C Storage, Dynamic Range 400 IU/mL - 4,000,000 IU/mL	IVD
CMVP200 CMV Verification Panel 1.4 mL	6 Panel Members, 1.4 mL, -20 °C Storage Dynamic Range 200 IU/mL - 4,000,000 IU/mL	IVD
CMVL101 CMV Low Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	CE-IVD, IVD
CMVH102 CMV High Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	CE-IVD, IVD
CMVB200 CMV Verification Bundle	4 CMVP200 Verification Panels, 1 Kit 1E3 IU/mL Precision (15 Vials), 1 Kit 5E5 IU/mL Precision (15 Vials), 2 Kits CMV LOD (10 Vials)	IVD
EBVP100 EBV Verification Panel	5 Panel Members, 1.0 mL, -20°C Storage, Dynamic Range 500 IU/mL - 5,000,000 IU/mL	CE-IVD, IVD
EBVL101 EBV Low Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	CE-IVD, IVD
EBVH102 EBV High Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	CE-IVD, IVD
HV6AP100 HHV-6A Verification Panel	5 Panel Members, 1.0 mL, -20°C Storage, Dynamic Range 500 cp/mL - 5,000,000 cp/mL	CE-IVD, IVD
HV6BP100 HHV-6B Verification Panel	5 Panel Members, 1.0 mL, -20°C Storage, Dynamic Range 500 cp/mL - 5,000,000 cp/mL	CE-IVD, IVD

Please note that not all products may be available in all countries.

SKU/PRODUCT NAME	DESCRIPTION	INTENDED USE
TRANSPLANTATION		
HV6L101 HHV-6 Low Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	CE-IVD, IVD
HV6H102 HHV-6 High Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	CE-IVD, IVD
HSV1P100 HSV-1 Verification Panel	5 Panel Members, 1.0 mL, -20°C Storage, Dynamic Range 200 cp/mL - 2,000,000 cp/mL	CE-IVD, IVD
HSV1L101 HSV-1 Low Run Control	5 Vials, 1.0 mL, Synthetic CSF Matrix, -20°C Storage	CE-IVD, IVD
HSV1H102 HSV-1 High Run Control	5 Vials, 1.0 mL, Synthetic CSF Matrix, -20°C Storage	CE-IVD, IVD
HSV2P100 HSV-2 Verification Panel	5 Panel Members, 1.0 mL, -20°C Storage, Dynamic Range 200 cp/mL - 2,000,000 cp/mL	CE-IVD, IVD
HSV2L101 HSV-2 Low Run Control	5 Vials, 1.0 mL, Synthetic CSF Matrix, -20°C Storage	CE-IVD, IVD
HSV2H102 HSV-2 High Run Control	5 Vials, 1.0 mL, Synthetic CSF Matrix, -20°C Storage	CE-IVD, IVD
HSVB200 HSV Verification Bundl	5 HSVP200 Verification Panels (9 Vials), 1.0 mL, -20°C Storage	IVD
JCVP100 JCV Verification Panel	6 Panel Members, 1.0 mL, -20°C Storage, Dynamic Range 200 IU/mL - 20,000,000 IU/m	CE-IVD, IVD
JCVL101 JCV Low Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	CE-IVD, IVD
JCVH102 JCV High Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	CE-IVD, IVD
VZVP100 VZV Verification Panel	5 Panel Members, 1.0 mL, -20°C Storage, Dynamic Range 500 cp/mL - 2,000,000 cp/mL	CE-IVD, IVD
VZVL101 VZV Low Run Control	5 Vials, 1.0 mL, Synthetic CSF Matrix, -20°C Storage	CE-IVD, IVD
VZVH102 VZV High Run Control	5 Vials, 1.0 mL, Synthetic CSF Matrix, -20°C Storage	CE-IVD, IVD



SKU/PRODUCT NAME	DESCRIPTION	INTENDED USE
VIROLOGY		
HBVP100 HBV Verification Panel 1.0 mL	7 Panel Members, 1.0 mL, -20°C Storage, Dynamic Range 100 IU/mL - 100,000,000 IU/mL	IVD
HBVP200 HBV Verification Panel 1.4 mL	7 Panel Members, 1.4 mL, -20°C Storage, Dynamic Range 100 IU/mL - 100,000,000 IU/mL	IVD
HBVL101 HBV Low Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	IVD
HBVH102 HBV High Run Contr	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	IVD
HBVB200 HBV Verification Bundle	4 HBVP200 Verification Panels, 1 Kit 1E3 IU/mL Precision (15 Vials), 1 Kit 1E6 IU/mL Precision (15 Vials), 2 Kits HBV LOD (10 Vials)	IVD
HCVP100 HCV Verification Panel 1.0 mL	7 Panel Members, 1.0 mL, -20°C Storage, Dynamic Range 100 IU/mL - 30,000,000 IU/mL	IVD
HCVP200 HCV Verification Panel 1.4 mL	7 Panel Members, 1.4 mL, -20°C Storage, Dynamic Range 100 IU/mL - 30,000,000 IU/mL	IVD
HCVL101 HCV Low Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	IVD
HCVH102 HCV High Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	IVD
HCVB200 HCV Verification Bundle	4 HCVP200 Verification Panels, 1 Kit 1E3 IU/mL Precision (15 Vials), 1 Kit 1E6 IU/mL Precision (15 Vials), 2 Kits HCV LOD (10 Vials)	IVD
HCVGTP HCV Genotype Panel	GT1a: 20,000 IU/mL, GT1b: 20,000 IU/mL, GT2: 20,000 IU/mL, GT3: 20,000 IU/mL, GT4: 20,000 IU/mL, GT5: 20,000 IU/mL	RUO
HIVP100 HIV-1 Verification Panel 1.0 mL	7 Panel Members, 1.0 mL, -20°C Storage, Dynamic Range 100 cp/mL - 6,000,000 cp/mL	IVD
HIVP200 HIV-1 Verification Panel 1.4 mL	7 Panel Members, 1.4 mL, -20°C Storage, Dynamic Range 100 cp/mL - 6,000,000 cp/mL	IVD
HIVL101 HIV-1 Low Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	IVD
HIVH102 HIV-1 High Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	IVD
HIVB200 HIV-1 Verification Bundle	HIVP200 Verification Panels, 1 Kit 1E3 cp/mL Precision (15 Vials), 1 Kit 1E5 cp/mL Precision (15 Vials), 2 Kits HIV-1 LOD (10 Vials)	IVD
AMRP100 Analytical Measurement Range Bundle	1 Panel of each HBVAMR, HCVAMR, HIVAMR	IVD
HBVAMR HBV Analytical Measurement Range Panel	3 Panel Members, 100 IU/mL, 100,000 IU/mL, 10,000,00 IU/mL, 1.0 mL, -20°C Storage	IVD
HCVAMR HCV Analytical Measurement Range Panel	3 Panel Members, 500 IU/mL, 50,000 IU/mL, 5,000,000 IU/mL, 1.0 mL, -20°C Storage	IVD
HIVAMR HIV-1 Analytical Measurement Range Panel	3 Panel Members, 600 cp/mL, 60,000 cp/mL, 6,000,000 cp/mL, 1.0 mL, -20°C Storage	IVD

Please note that not all products may be available in all countries.

SKU/PRODUCT NAME	DESCRIPTION	INTENDED USE
VIROLOGY		
ZKVP100 ZKV Verification Panel	5 Panel Members, 1.0 mL, -70°C Storage, Dynamic Range 100 cp/mL - 1,000,000 cp/mL	CE-IVD, IVD
ZKVL101 ZKV Low Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -70°C Storage	CE-IVD, IVD
ZKVH102 ZKV High Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -70°C Storage	CE-IVD, IVD
NEGATIVE		
NEG000 Negative Run Control	5 Vials, 1.0 mL, True EDTA Plasma, -20°C Storage	IVD
NEG050 Bulk Negative	50 mL Bottle, True EDTA Plasma, -20°C Storage	IVD
HPV		
HPVP100 HPV Control Set (HPV GT 16, 18, 68, and Negative)	1 Kit of each: HPV16C, HPV18C, HPV68C, HPVNEG, 10 Vials per kit, 5.0 mL, PreservCyt, 2-8°C Storage	CE-IVD, IVD
HPV16C HPV Genotype 16 Control	10 Vials, 5.0 mL, PreservCyt, 2-8°C Storage	CE-IVD, IVD
HPV18C HPV Genotype 18 Control	10 Vials, 5.0 mL, PreservCyt, 2-8°C Storage	CE-IVD, IVD
HPV68C HPV Genotype 68 Control	10 Vials, 5.0 mL, PreservCyt, 2-8°C Storage	CE-IVD, IVD
HPVNEG HPV Negative Control	10 Vials, 5.0 mL, PreservCyt, 2-8°C Storage	CE-IVD, IVD
MICROBIOLOGY		
CDFB200 C. difficile Verification Bundle	5 CDFP200 Verification Panels (3 Vials), 1.0 mL, -20°C Storage	CE-IVD, IVD
CTNGRC CT/NG Run Control	6.0 mL Dropper, Synthetic Matrix, 2-8°C Storage	IVD
CTNGB200 CT/NG Verification Bundle	5 CTNGP200 Verification Panels (9 Vials), 1.0 mL, -20°C Storage	IVD
MRSAB200 MRSA Verification Bundle	5 MRSAP200 Verification Panels (6 Vials), 1.0 mL, -20°C Storage	CE-IVD, IVD



SKU/PRODUCT NAME	DESCRIPTION	INTENDED USE
MENINGITIS		
EVRC Enterovirus Run Control	5 Vials, 1.0 mL, Synthetic Matrix, -20°C Storage	CE-IVD, IVD
NEW / RESPIRATORY		
RPRC Respiratory Run Control	2 Vials of Each: Mix 1, Mix 2, Mix 3, Mix 4 - 0.3 mL Fill Volume, -20°C Storage	CE-IVD, IVD
RRPOS Respiratory Positive Run Control	6 Vials, 0.3 mL Fill Volume, -20°C Storage	CE-IVD, IVD
RPNEG Respiratory Negative Run Control	6 Vials, 0.3 mL Fill Volume, -20°C Storage	CE-IVD, IVD
COV019 SARS-CoV-2 Standard	5 Vials, 0.3 mL Fill Volume, -20°C or below Storage, E, N and S genes of SARS-CoV-2, 200,000 cp/mL each	RUO
COV000 SARS-CoV-2 Negative	5 Vials, 0.3 mL Fill Volume, -20°C or below Storage 75,000 cp/mL	RUO
COV019CE SARS-CoV-2 Positive Run Control	5 Vials, 0.3 mL Fill Volume, -20°C or below Storage, E, N and S genes of SARS-CoV-2, 200,000 cp/mL each	CE-IVD, IVD
COV000CE SARS-CoV-2 Negative Run Control	5 Vials, 0.3 mL Fill Volume, -20°C or below Storage 75,000 cp/mL	CE-IVD, IVD
VECTOR BORNE		
ANARC Anaplasma Run Control	5 Vials, 1.0 mL, Synthetic Matrix, -20° Storage	CE-IVD, IVD
BABRC Babesia Run Control	5 Vials, 1.0 mL, Synthetic Matrix, -20° Storage	CE-IVD, IVD
BRTRC Bartonella Run Control	5 Vials, 1.0 mL, Synthetic Matrix, -20° Storage	CE-IVD, IVD
BORRC Borrelia Run Control	5 Vials, 1.0 mL, Synthetic Matrix, -20° Storage	CE-IVD, IVD
ERLRC Ehrlichia Run Control	5 Vials, 1.0 mL, Synthetic Matrix, -20° Storage	CE-IVD, IVD

Please note that not all products may be available in all countries.

Custom Manufacturing



EXACT Diagnostics Development Stage Features

Discovery & Feasibility

- Infectious Disease, Genetics, and Oncology materials
- Esoteric analytes and matrices available
- Value assignment utilizing Real Time and Digital PCR
- International Standards available

IVD Development

- Panels to assess your dynamic range, analytical sensitivity and specificity
- Manufacture and prepare specimens for clinical trials

Manufacturing

- All Products manufactured under cGMP
- Customizable OEM standards and control materials
- Quality Control materials for batch release

510(k) / PMA

- Data Analysis / Regulatory Support

Market / Commercialization

- Catalog products assist manufacturer customers in validation and daily run controls
- Operator and Probit Analysis Panels
- IVD materials for your IVD Assay



A Bio-Rad Laboratories Company

For more information or ordering, contact:
CustomerService@exactdiagnostics.com or call 817.989.9262

Exact Diagnostics, 3400 Camp Bowie Blvd. CBH-214, Fort Worth, Texas 76107.