

Ensure sensitive host cell contaminant detection

Be safe in quantifying residual host cell DNA and proteins when developing your next biologic, cell or gene therapy

CHO, HEK293, and E. coli cells are common hosts for bioproduction of antibody-based drugs and cell and gene therapies (CGT). Product impurities from residual host cell proteins (HCP) or host cell DNA need to be minimized, as too high levels raise safety concerns due to potential immunogenic and oncogenic effects.

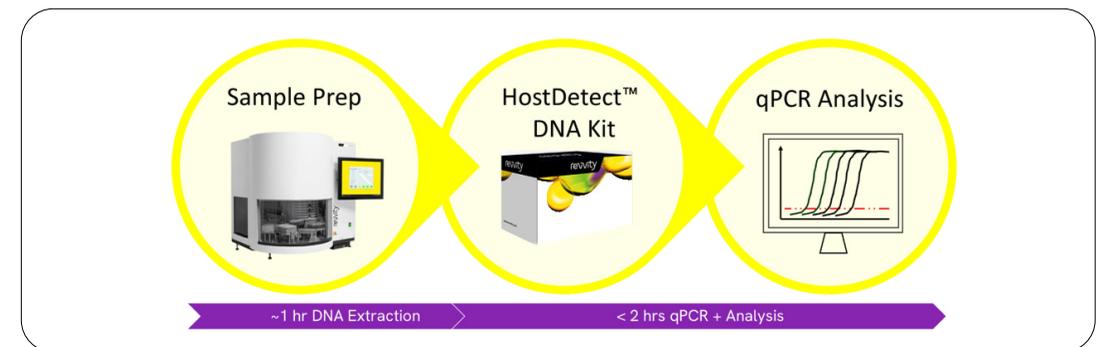
Elevate safety standards with residual DNA detection solutions HostDetect™

Our bioprocess quality control solutions are designed to streamline host cell residual DNA detection workflow, from sample extraction to PCR quantification, to ensure the purity of biotherapeutic or CGT products via demonstrating that host cell contaminant levels are in compliance with critical to quality attributes.

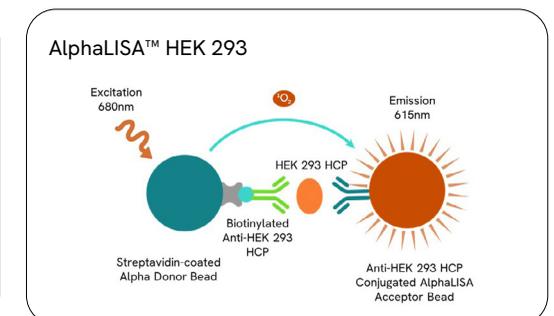
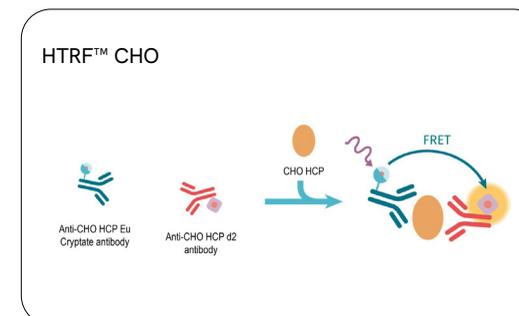
Monitor host cell protein impurities with HTRF™ and AlphaLISA™

Our HTRF and AlphaLISA CHO & HEK293 HCP assays enable ready-to-use homogeneous, no-wash detection and quantification of CHO and HEK293 HCP impurities. These off-the-shelf kits deliver a streamlined workflow, a broader dynamic range and higher analytical sensitivity than traditional multi-step ELISA assays.

DNA detection



HCP detection



HostDetect™ streamlines your workflows of residual DNA testing of biologics

Our HostDetect CHO (DXMDX-RGT-1003), HEK293 (DXMDX-RGT-1004), and E.coli (DXMDX-RGT-1005) PCR DNA Quant Kits integrates chemagic™ DNA extraction and real-time PCR technologies to provide a seamless workflow from sample to result in less than 3 hours. Benefit from a streamlined detection process that helps to accelerate biologics or CGT product development and purification timelines, as well as enhancing purity and safety.

Precise contaminant high throughput DNA quantification via automation

- Up to 96 samples from in-process or purified final products with different matrices.
- Consistent performance and increased sample processing throughput

Detect and quantify residual DNA with increased precision and analytical sensitivity

- Utilize TaqMan-probe based qPCR with a qualified DNA reference standard.
- Internal control provided to verify experiment procedures and assay integrity
- Documentation and support includes limit of quantitation (LOQ), analytical specificity and accuracy, continuous updates to align with evolving residual DNA testing regulations

Customizing tailored solutions

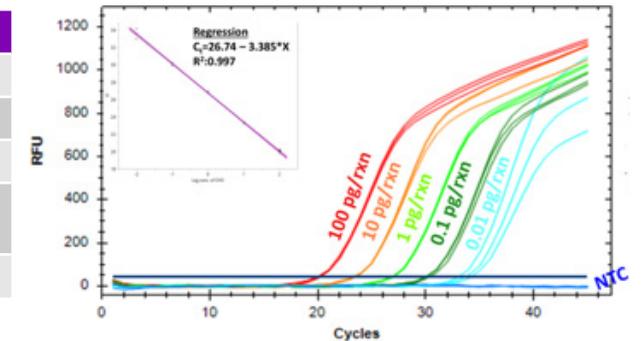
- Utilize your own established nucleic acid extraction techniques for sample preparation and operate a range of conventional qPCR devices.
- Comprehensive technical expertise for tailored solutions to meet your specific requirements

HTRF and AlphaLISA kits for reliable host cell protein impurity quantification

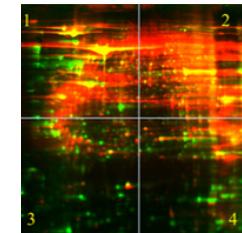
Our no-wash HTRF (64CHOPEG/H) and AlphaLISA (AL3198HV/C/F and AL3176HV/C/F) HCP kits quantify residual CHO and HEK293 proteins with enhanced precision, analytical sensitivity, and efficacy.

- User-friendly with no-wash steps
- Easily automatable
- Fewer touchpoints for reproducible handling
- Broad dynamic range
- Heightened analytical sensitivity
- Assays display exceptional dilutional linearity
- Compatible with the most used biologics manufacturing buffers
- Reliable antigen spike recovery
- Robust reproducibility
- Excellent HCP coverage

	HostDetect CHO
LOQ	0.01 pg/mL (5pg/mL sample)
Range	100 pg to 0.01 pg/rxn
Linearity	$R^2 \geq 0.98$
Analytical specificity	Negative of HEK293, E.coli
Precision	CV < 5% at 0.01 pg/rxn



	AlphaLISA - HEK293	AlphaLISA - CHO
LOQ (ng/mL)	1.01	1.80
Range (ng/mL)	1.01-1000	1.80-3,000
Precision	CV < 15%	
Time to results	3h30	



HEK 293 HCPs (green) coverage of anti-HEK293 (red) polyclonal antibody = 98.7%

Learn more about our residual DNA and HCP detection and quantification solutions.

revvity