

EnVision® Nexus™ multimode plate reader

The EnVision Nexus comes with filter-based technology for absorbance, fluorescence, luminescence, TRF, and fluorescence polarization. We've combined filters and optics into a single module optimized for your assays. These new filter modules are held in exact position magnetically and can be exchanged without tools, making them simple to install and replace. We've integrated these filter modules with Kaleido, so as soon as you load a new module, the software recognizes it.

To find the filter module optimized for your applications, please search the list below by measurement technology, wavelength, or assay to see the part number we recommend. Compatible fluorophores are listed next to each filter module where relevant.

Note: This is not an exhaustive list of relevant fluorophores. Please reach out to your local Revvity representative for support when selecting the right filter module for your application.

Fluorescence Intensity

Description	Part Number	Channel	Excitation Filter	Emission Filter (1 + 2)	Dichroic Mirror (1 + 2)	Additional Fluorophores
FITC PicoGreen	HH36791001	1	485	535	D509	Alexa 488, PhenoVue488, Calcein, DyLight 488, FAM, Fluo-4, (e)GFP, Rhodamine 110, SybrGreen
Beta Lactamase	HH36791002	2	400	535 + 460	D425 + D490	
Phenovue 647	HH36791003	1	632	685	D658	Alexa 647, Cy5
Phenovue 568	HH36791004	1	535	615	D585	Alexa 568, Texas Red
Hoechst	HH36791005	1	355	450	D409	PhenoVue Hoechst 33342
Umbelliferone	HH36791007	1	355	460	D409	Alexa 350, AMC , BFP, DAPI
Texas Red	HH36791008	1	555	632	D595	mCherry, ROX, PhenoVue 594, Alexa 594
Bodipy	HH36791011	1	535	595	D555	Amplex Red, Resorufin
Alamar Blue	HH36791012	1	530	580	D555	PhenoVue555, 5-TAMRA, Cy3, HEX, mOrgange, Td Tomato, TRITC
Fura-2	HH36791013	1	340	510	D409	

Fluorescence Polarization

Description	Part Number	Channel	Excitation Filter	Emission Filter (1 + 2)	Dichroic Mirror (1 + 2)	Additional Fluorophores
FP FITC	HH36792001	2	485	535p + 535s	D509 + D535fp	Alexa 488, PhenoVue488, Calcein, DyLight 488, FAM, Fluo-4, (e)GFP, Rhodamine 110, SybrGreen
FP Cy5	HH36792002	2	620	685p + 685s	D658 + D685fp	Alexa 647, APC, Phenovue 647
Texas Red FP	HH36792004	2	555	632p + 632s	D595 + D632fp	Alexa 568, Alexa 594 PhenoVue 594
Bodipy FP	HH36792005	2	535	595p + 595s	D555 + D595fp	

Luminescence

Description	Part Number	Channel	Excitation Filter	Emission Filter (1 + 2)	Dichroic Mirror (1 + 2)
General Lum	HH36793001	1	N/A	700	N/A
NanoBRET	HH36793002	2	N/A	645 + 460	D585
BRET	HH36793003	2	N/A	530 + 485	D509
BRET2	HH36793004	2	N/A	515 + 410	D475
ChromaGlo	HH36793005	2	N/A	645 + 510	D585

Time Resolved Fluorescence (TRF)

Description	Part Number	Channel	Excitation Filter	Emission Filter (1 + 2)	Dichroic Mirror (1 + 2)
DELFI A	HH36794001	1	330	615	D409
HTRF LANCE	HH36794002	2	330	665 + 620	D409 + D645
HTRF LANCE Las.	HH36794003	2	N/A	665 + 620	D409 + D645
HTRF Green	HH36794004	2	330	620 + 515	D409 + D555
HTRF Green Las.	HH36794005	2	N/A	620 + 515	D409 + D555
DELFI A Eu Las.	HH36794006	1	N/A	615	D409

Alpha

Description	Part Number	Channel	Excitation Filter	Emission Filter (1 + 2)	Dichroic Mirror (1 + 2)
Alpha Screen	HH36795001	1	N/A	575	D660
AlphaplexEu/Tb	HH36795002	2	N/A	615 + 546	D660 + D585
Alphaplex Eu/Sm	HH36795003	2	N/A	644 + 615	D660 + D630
Alphaplex Sm/Tb	HH36795004	2	N/A	644 + 546	D660 + D585

Absorbance

Description	Part Number	Channel	Excitation Filter	Emission Filter (1 + 2)	Dichroic Mirror (1 + 2)
ELISA ABTS	HH36796001	1	405	N/A	Mirror
ELISA TMB	HH36796002	1	450	N/A	Mirror
Nucl. Ac. Abs.	HH36796005	1	260	N/A	Mirror
Prot. Abs.	HH36796006	1	280	N/A	Mirror
ELISA OPD	HH36796007	1	492	N/A	Mirror
Quinoneimine dy	HH36796008	1	540	N/A	Mirror
BCA Prot. assay	HH36796009	1	560	N/A	Mirror
MTT Assay	HH36796010	1	570	N/A	Mirror
Bradford assay	HH36796011	1	595	N/A	Mirror
Bacteria Abs.	HH36796012	1	600	N/A	Mirror
Lowry Prot. Ass	HH36796013	1	750	N/A	Mirror
Pierce Prot Ass	HH36796014	1	660	N/A	Mirror

