

NUNC™

Colour Coded Modules



Your Partner in
Immunology

Thermo
SCIENTIFIC

Contact us for Sales and Service
thermoscientific.com/contactus

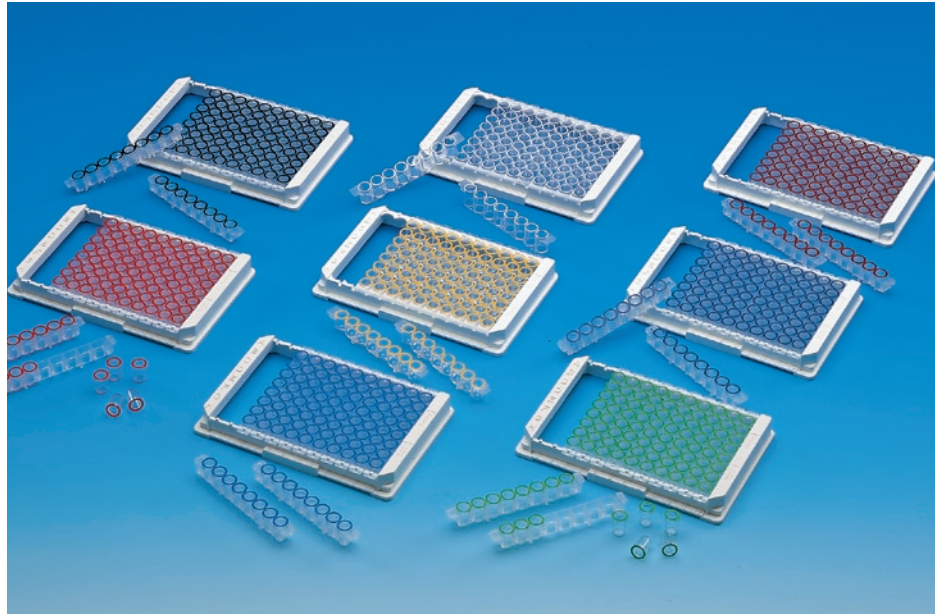
*Contact information contained within this document may be incorrect.

Nunc Colour Coded Modules

Nunc-Immuno Modules with colour coded rims are standardized to facilitate easy recognition of your assays and high reliability when working with individual wells.

This is a post-moulding process whereby a layer of pigment is applied to the plastic rim by heat stamping. Only one colour per plate is available. After the colour coding procedure the surface is retested in our laboratories to ensure that the binding properties conform to the required standard. All colour coated products include a certificate stating conformaty.

Before colour coding pre-batch approval is possible on request.



The colours are available in following formats:

BA = BreakApart, LW = LockWell

Colour	C8 MaxiSorp	C8 B/A MaxiSorp	C8 B/A PolySorp	C8 LW MaxiSorp	C8 LW PolySorp	U8 MaxiSorp	U8 PolySorp	F8 MaxiSorp
Ref. product.	445101	473768	473539	446469	446442	475078	475086	468667
Green	445120	459048	465005	475000	475443	478178	478191	462456
Yellow	445113	463630	464912	473237	475444	478176	478192	459990
Blue	445110	460798	460799	474691	475001	478179	478193	462455
Red	445117	460801	464904	475426	475445	475540	478194	463975
Dark Blue	445121	463649	465006	476422	475446	478180	478195	465560
White	445114	463657	465013	475442	475447	475541	478196	462465
Black	445118	465919	465007	475957	475454	478181	478202	462475
Orange	445115	465020	464920	475967	475448	478177	478197	465552
Burgundy	445119	459049	465008	477089	475449	478183	478198	462485
Grey	445112	463622	465009	478174	475451	478184	478199	462495
Brown	445111	465188	465010	478123	475452	478185	478200	462445
Purple	445116	465021	465011	475965	475453	478175	478201	468993
Turquoise	445122	465189	465012	476423	475455	478186	478205	462457
Maroon	445123	465190	465014	476424	475456	478187	478206	462458
Pink	445124	463469	465015	476425	475457	478188	478207	462459

Minimum order of two cases; all products are packed 10/60. Other modules can be colour coded - please ask for a quotation at your local Nunc distributor or at Infocity@nunc.dk.



Nunc A/S · Denmark · Tel +45 4631 2000 · infocity@nunc.dk · www.nuncbrand.com
 Nalge Nunc International · USA · Tel +1 585 586 8800 · nnitech@nalgennunc.com · www.nuncbrand.com
 Nunc GmbH & Co. KG · Germany · Tel +49 6111 86740 · nunc@nunc.de · www.nunc.de
 Nalge Nunc International KK · Japan · Tel +81 3381 63355 · info@nalgennunc.co.jp · www.nalgennunc.co.jp

www.nuncbrand.com