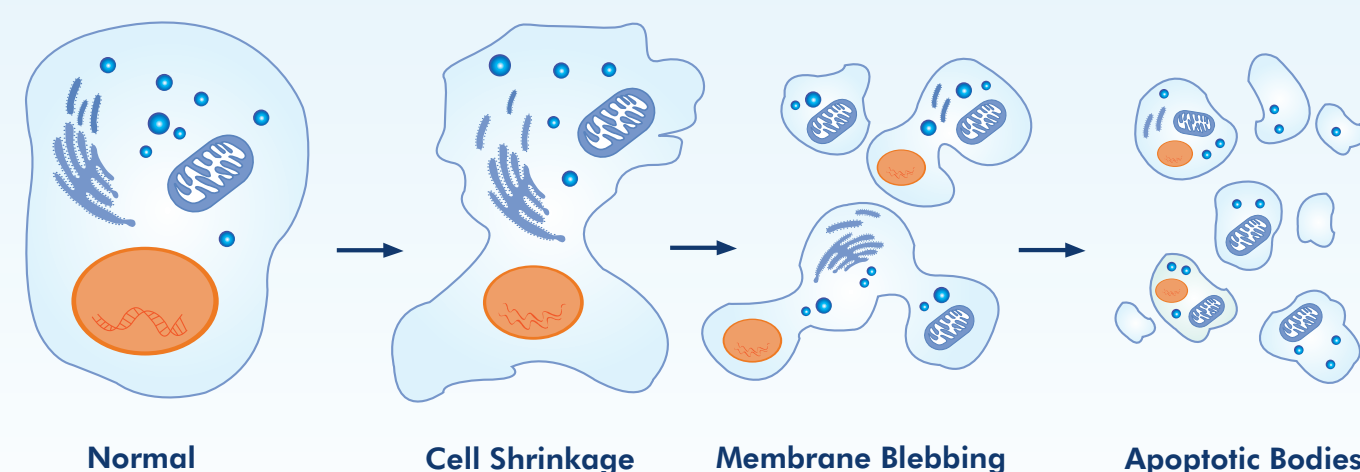
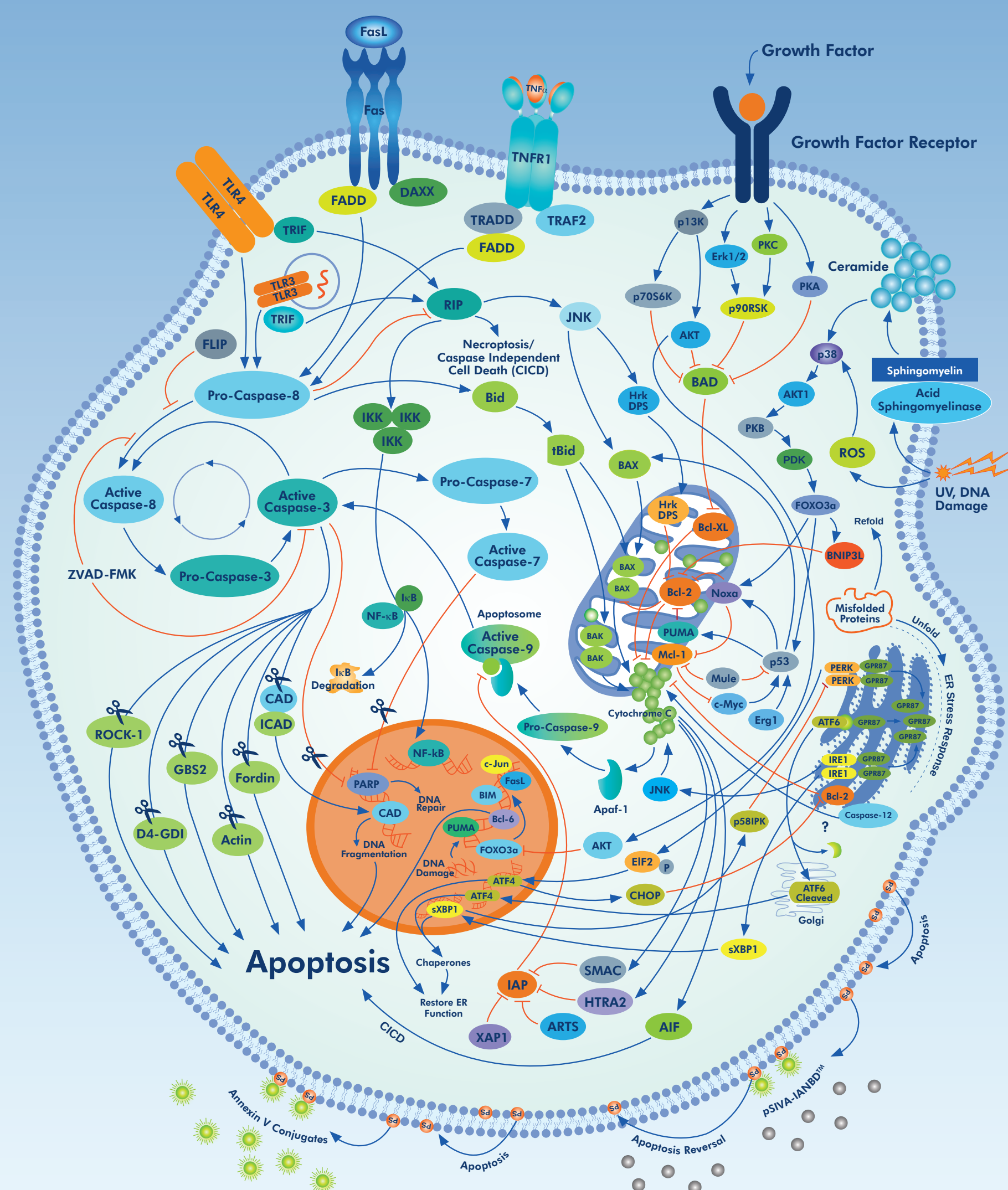
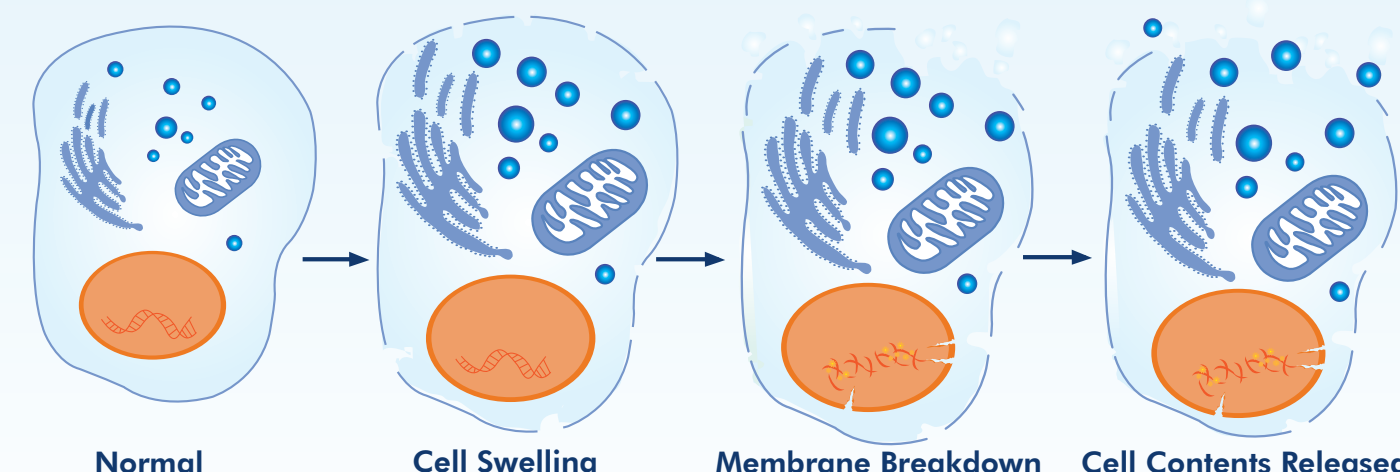
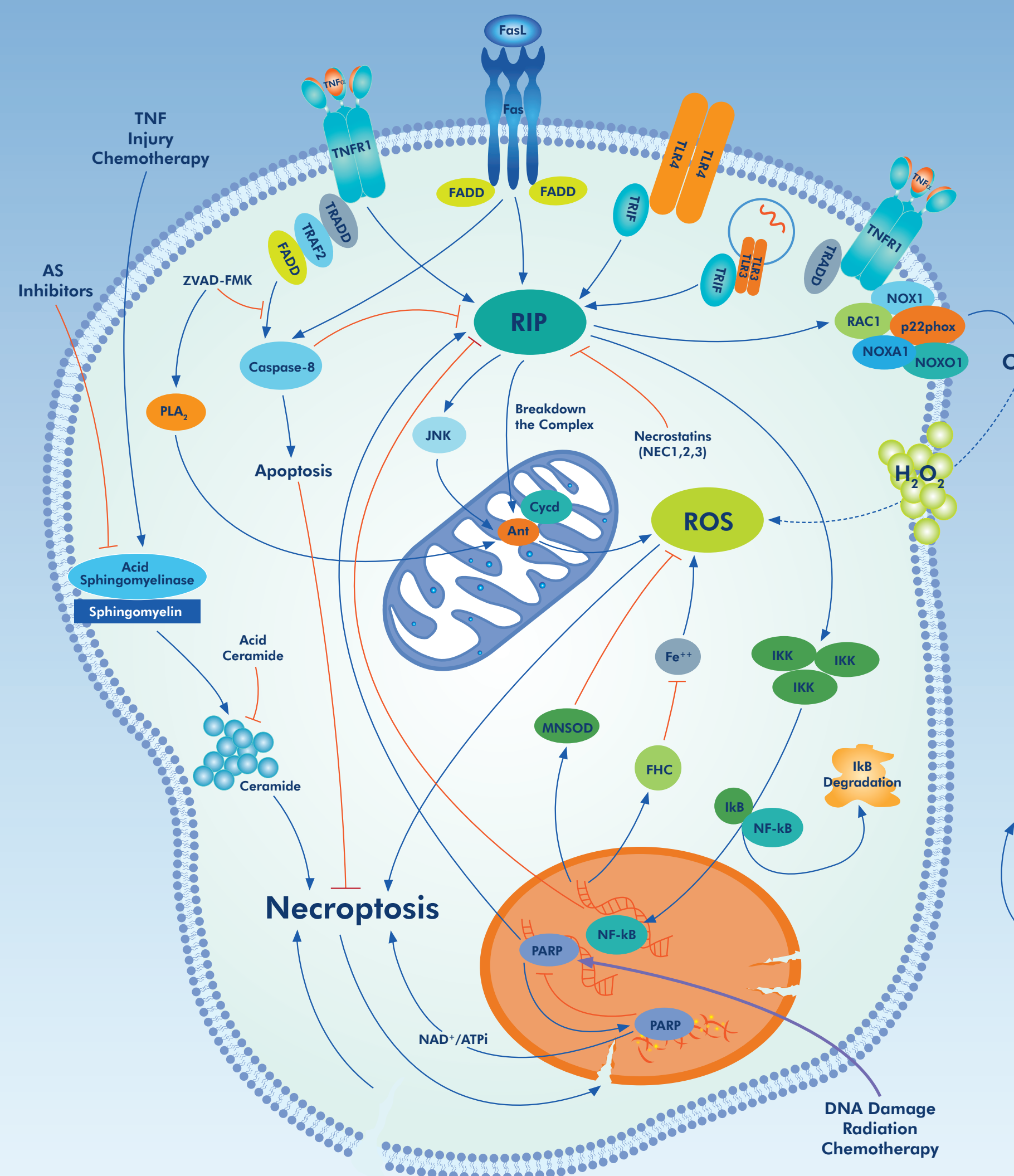


Apoptosis, Necroptosis & Autophagy

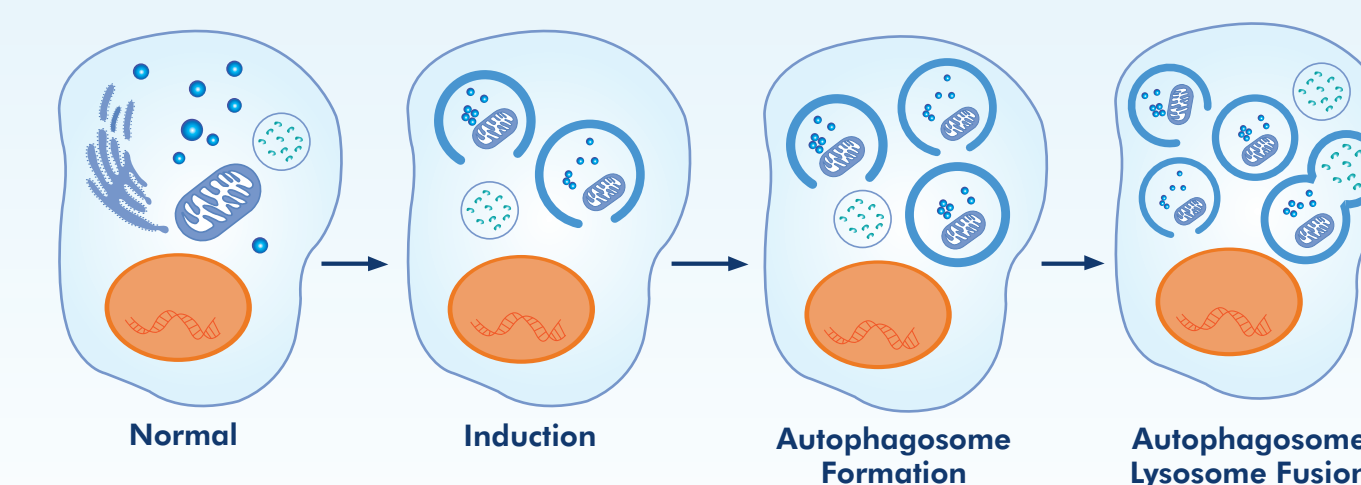
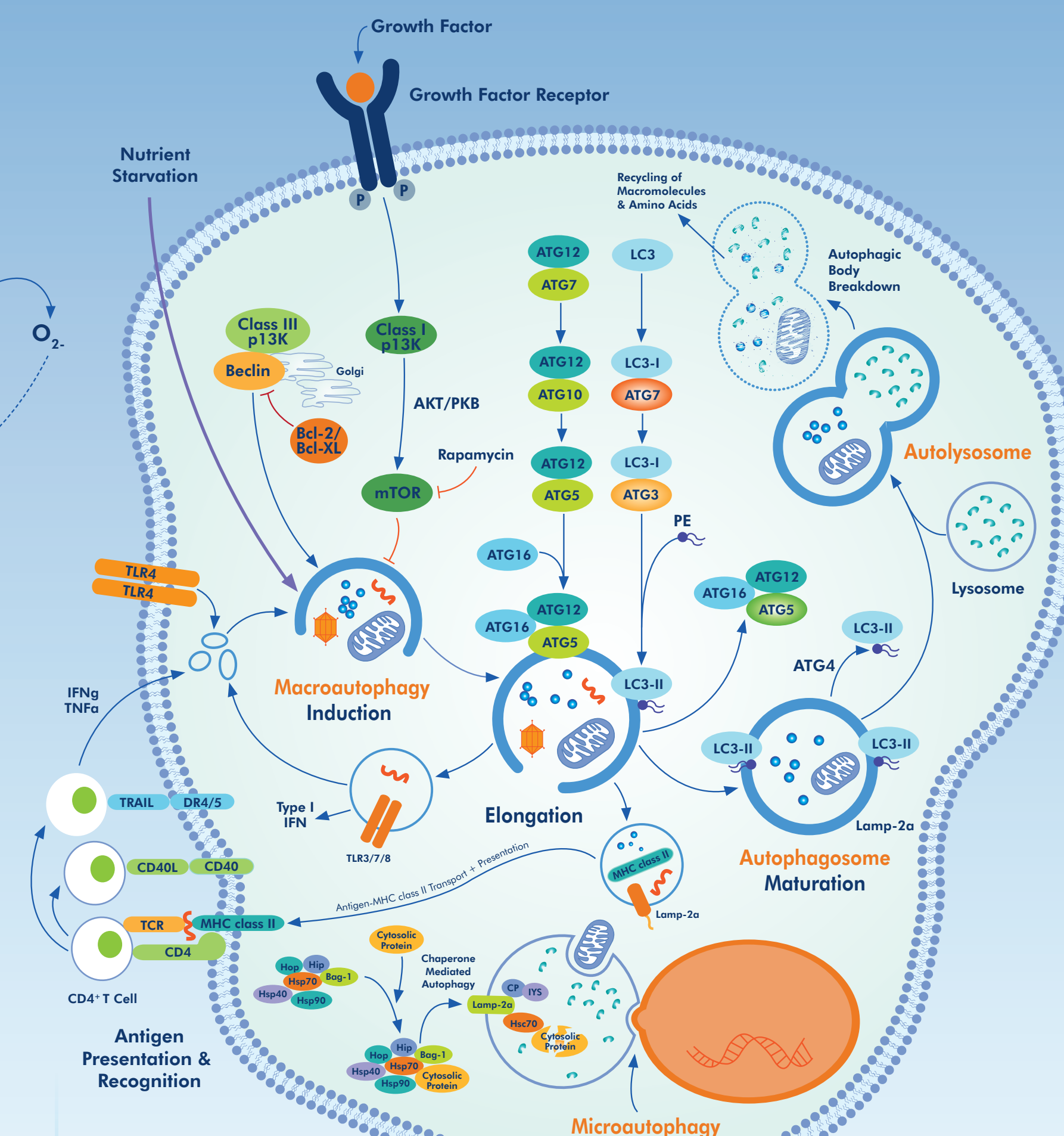
APOPTOSIS



NECROPTOSIS



AUTOPHAGY

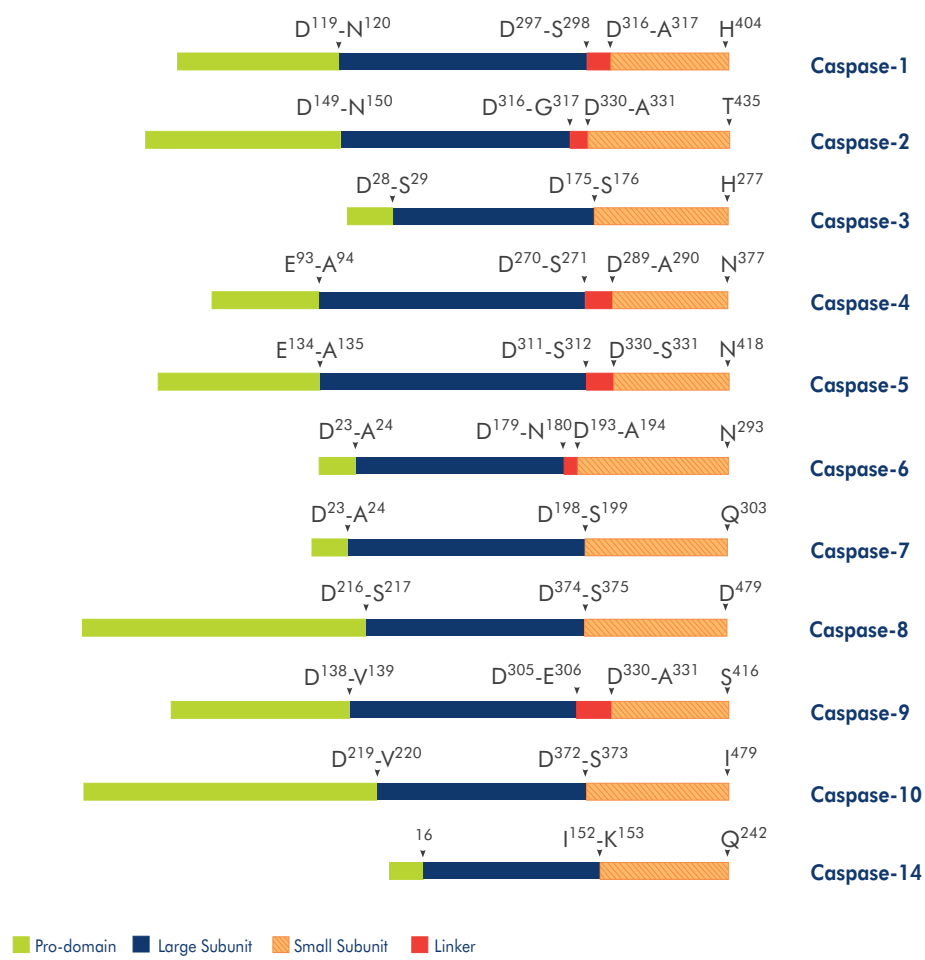


Cell Death Assays: According to the Recommendations of the Nomenclature Committee on Cell Death

The cell death field exploded during the 1990s and by the early 2000's assays to measure cell death were commonplace. New cell death types were routinely being elucidated and certain terms like 'apoptosis' became popularized. However descriptors like 'percent apoptosis' were vague, nomenclature lacked uniformity, and the cell death literature was becoming unwieldy. Hence in 2005, the editors of *Cell Death and Differentiation* formed the Nomenclature Committee on Cell Death (NCCD) which addressed the need to streamline cell death reporting. The NCCD guidelines are based on a series of measurable biochemical, functional and molecular features to characterize cell death subroutines. According to the NCCD, it is necessary to use at least two detection methods to precisely define an observed cell death.

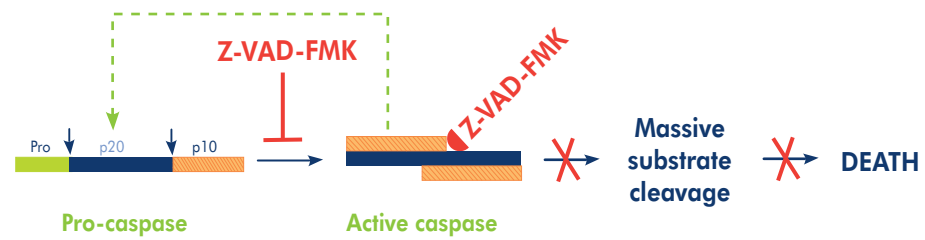
Novus Biologicals is proud to facilitate the recommendations of the NCCD by offering a broad, comprehensive portfolio encompassing a myriad of methods for studying cell death-related phenomenon.
Galluzzi et al. Cell Death Differ. 12:1463-1467 (2005). Kroemer et al. Cell Death Differ. 19:107-120 (2012).

Caspase Structures



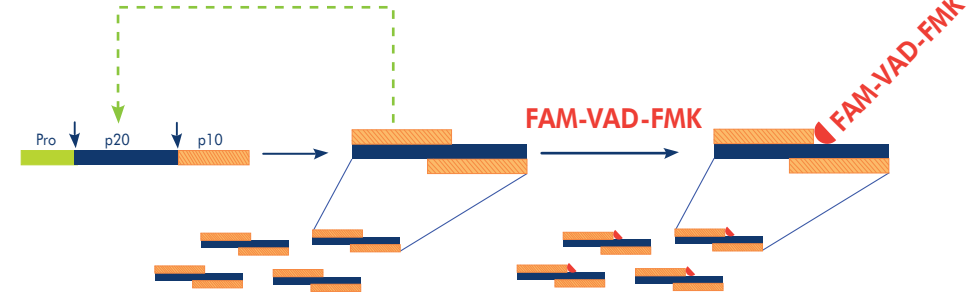
Caspases are synthesized as inactive pro-enzyme precursors containing an N-terminal prodomain of variable length followed large (p20) and small (p10) subunits. Caspases are activated through proteolytic cleavage (arrow symbols) at specific asparagine residues located within the pro-domain, and the p10 and p20 subunits. Activation results in the generation of mature active caspases consisting of heterotetramers p20₂p10₂.

Apoptosis Inhibition with Caspase Inhibitors



Caspase inhibitors like pan caspase inhibitor Z-VAD-FMK (NB22-29392) are used in vitro and in vivo assays to inhibit apoptosis. These cell permeable peptides function by binding irreversibly to the catalytic or active site of caspases, thereby preventing the processing of pro-caspases to their fully active forms.

Apoptosis Detection with Labeled Caspase Inhibitors



Fluorescently labeled caspase inhibitors like FAM-VAD-FMK (NB22-29383) are used in fluorescence microscopy and flow cytometry assays to detect apoptosis. Just like their unlabeled counterparts, they bind to the active site of caspases.

Caspase Inhibitors

Broad Spectrum: Boc-D(OMe)-FMK
Broad Spectrum: Z-VAD-(OMe)-FMK
Broad Spectrum: Q-VD-OPH
Caspase 3/7: Z-DEVD-FMK
Caspase 8: Z-IETD-FMK
Caspase 9: Z-LEHD-FMK
Negative Control: Z-FA-FMK

NB22-29395
NB22-29392
NB22-29391
NB22-29396
NB22-29397
NB22-29398
NB22-29384

Cell Death Terms

Accidental cell death
Anoikis
Aponecrosis
Apoptosis
Autolysis
Autophagic cell death
Autophagy

Caspase-dependent intrinsic apoptosis
Caspase-independent intrinsic apoptosis
Cornification
Entosis
Excitotoxicity
Extrinsic apoptosis by death receptors
Extrinsic apoptosis by dependence receptors

Mitotic catastrophe
Necrosis
Netosis
Necroapoptosis
Necroptosis
Parthanatos

Programmed cell death
Pyroptosis
Oncosis
Wallerian degeneration

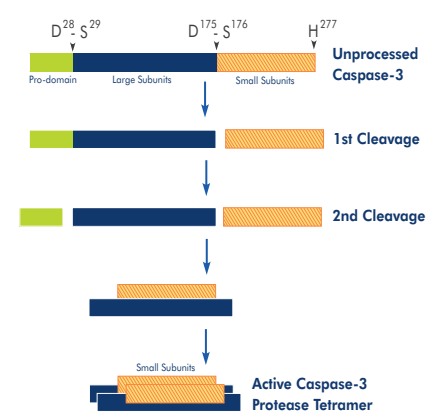
Antibody Detection of Pro and Active/Cleaved Caspases

Caspase antibodies are classical tools for detecting unprocessed (pro) and active (cleaved) forms of caspases by western blot. Western blot banding patterns will depend on the epitope recognized by a given antibody and the caspase forms present.

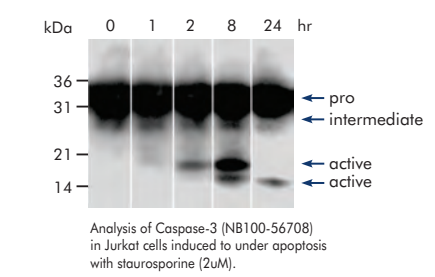
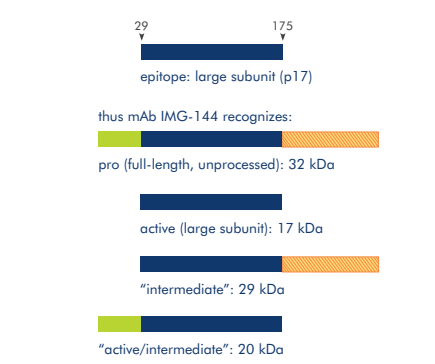
For example, the epitope of the Caspase-3 mAb (NB100-56708), clone 31A1067 has been mapped to the large subunit of Caspase-3 (between amino acids 29-176). Therefore, NB100-56708 recognizes pro-Caspase-3 (~32 kDa) and cleavage fragments containing the 17 kDa large subunit.

It is important to note that caspase cleavage fragments may be transient due to a short half-life and therefore loss of the caspase-3 pro-form without the appearance of cleavage fragments can also be an indication of caspase activation.

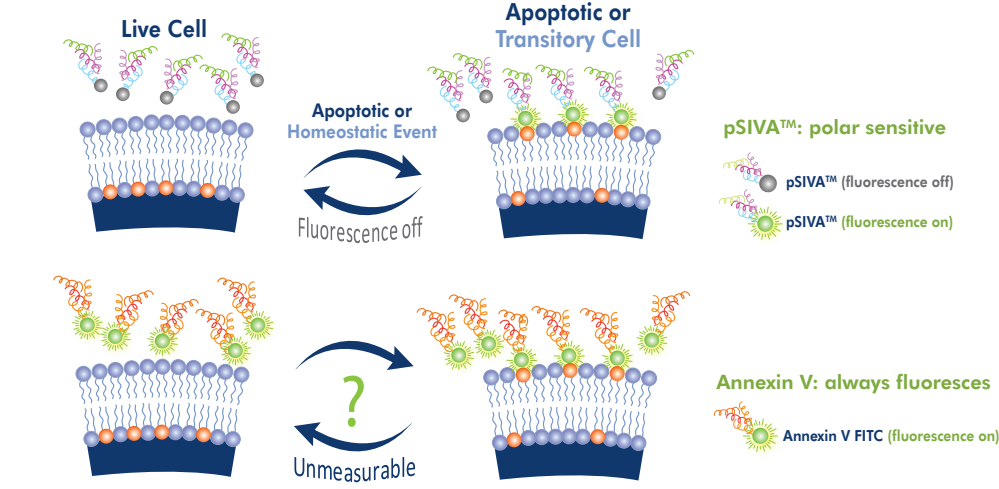
Caspase-3 Activation



Caspase-3 Antibody

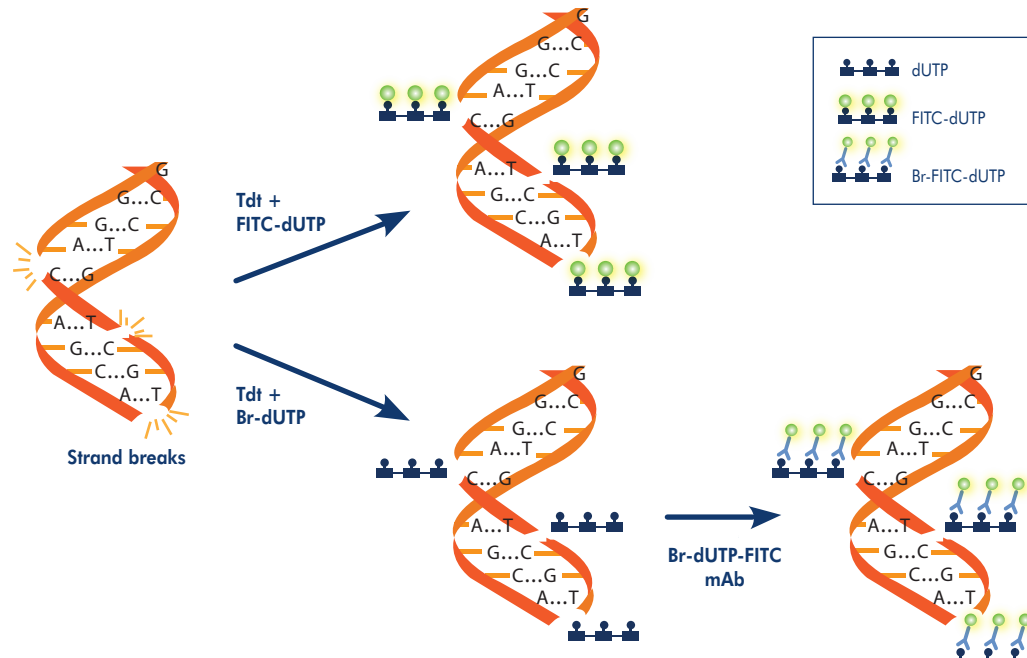


pSIVA-IANBD™ versus Annexin V-FITC



pSIVA-IANBD™ [Polarity Sensitive Indicator of Viability (NB22-29382/NB22-29611)] and Annexin V-FITC (NB22-29373) apoptosis technology. Both pSIVA-IANBD™ and Annexin V-FITC fluorescent conjugates bind to exposed bind to exposed phosphatidylserine (PS) residues on cell surfaces, a hallmark of apoptosis. A key advantage of pSIVA-IANBD™ is that it fluoresces only when bound to cells, rendering it uniquely suitable for live microscopy imaging of PS exposure and monitoring apoptosis real time.

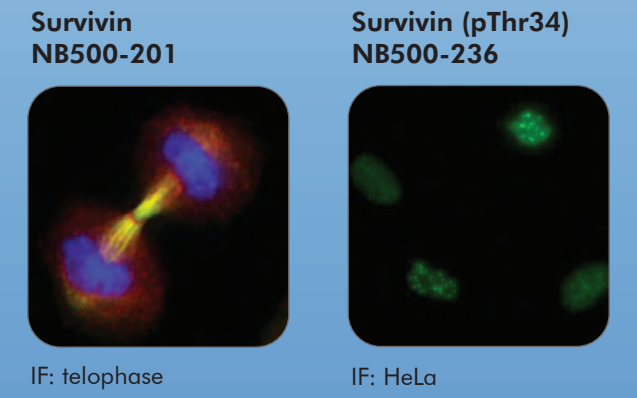
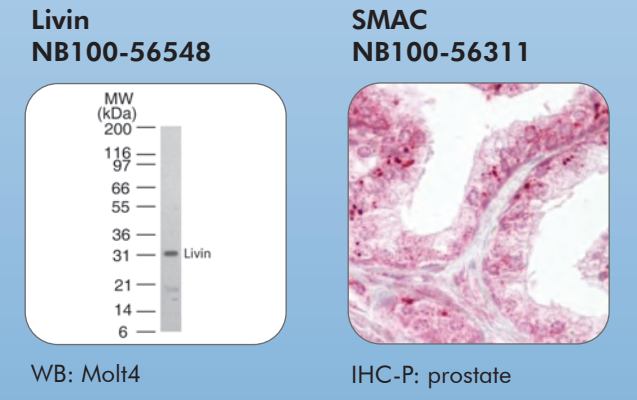
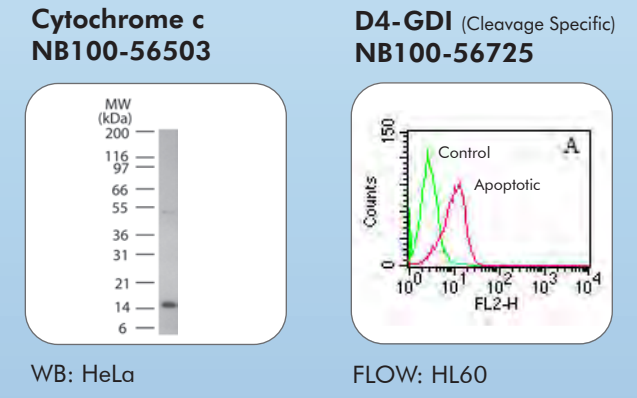
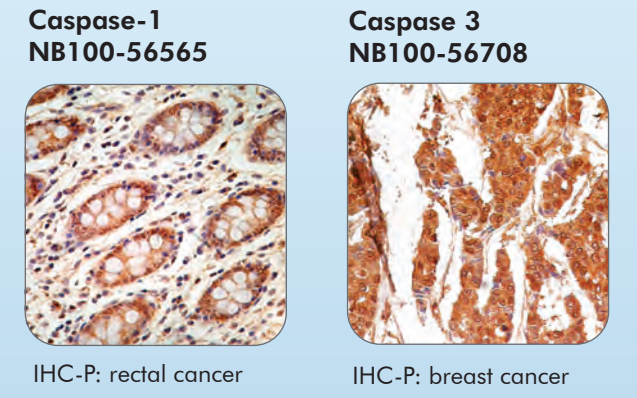
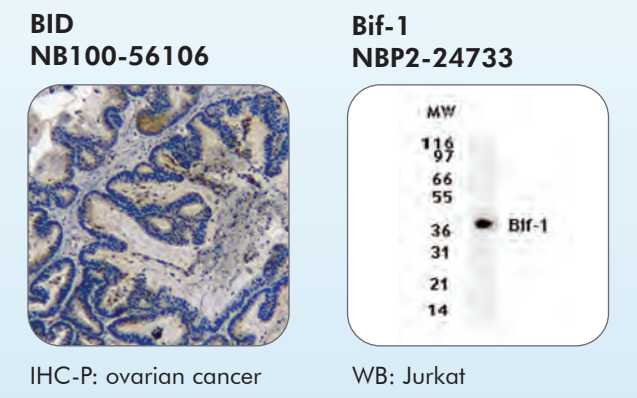
DNA Fragmentation TUNEL Apoptosis Assays



APO-BRDU™ (NB22-31161, NB22-31164) and APO-DIRECT™ (NB22-31159) TUNEL Assays detect DNA fragmentation associated with apoptosis. TdT catalyzes the addition of Br-dUTP or FITC-dUTP to DNA strand breaks. FITC-dUTP is detected directly and Br-dUTP is detected with a BrdU-FITC mAb.

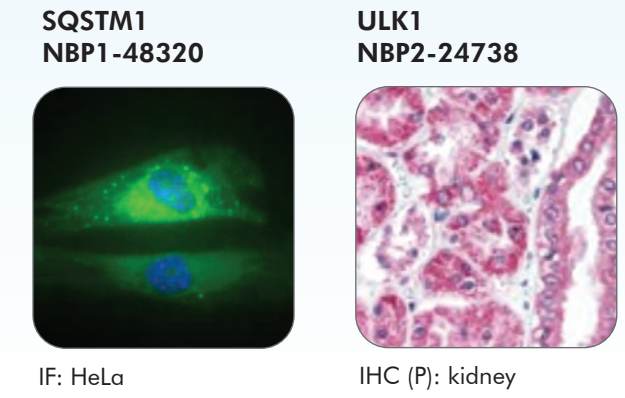
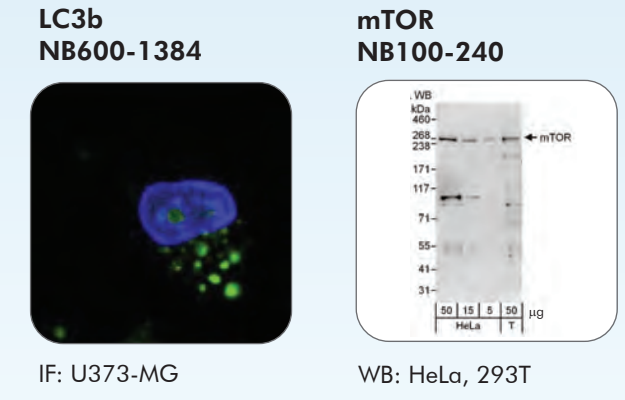
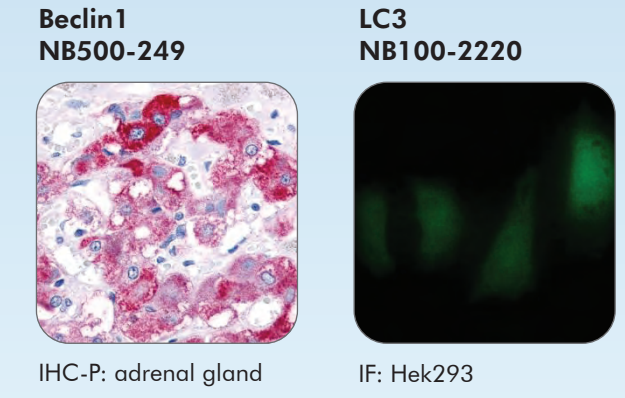
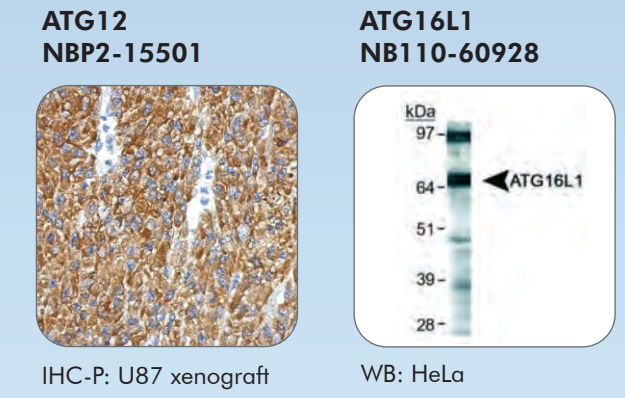
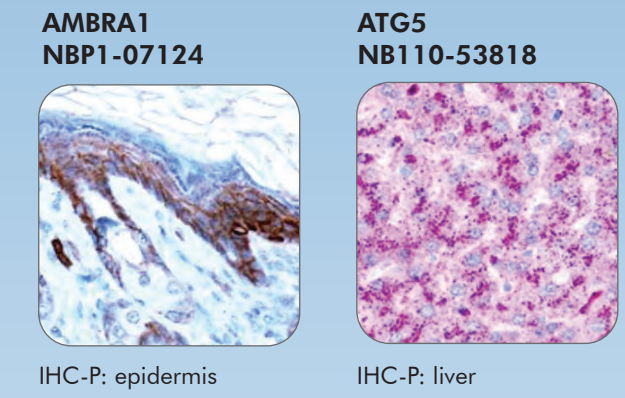
Apoptosis

Apoptosis is a naturally occurring, programmed cell death process involving a series of biochemical events leading to characteristic morphological changes and cellular demise. Apoptosis dysregulation has been implicated in a myriad of diseases.



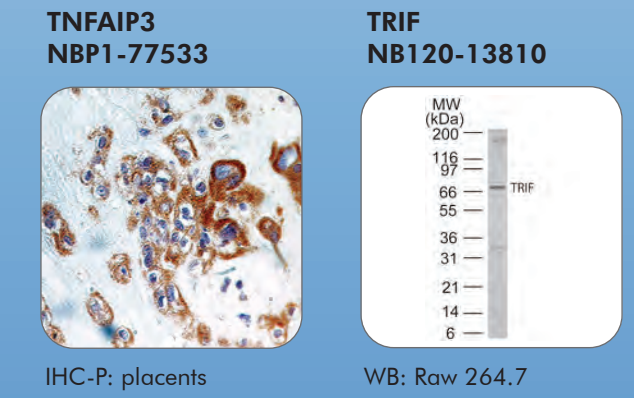
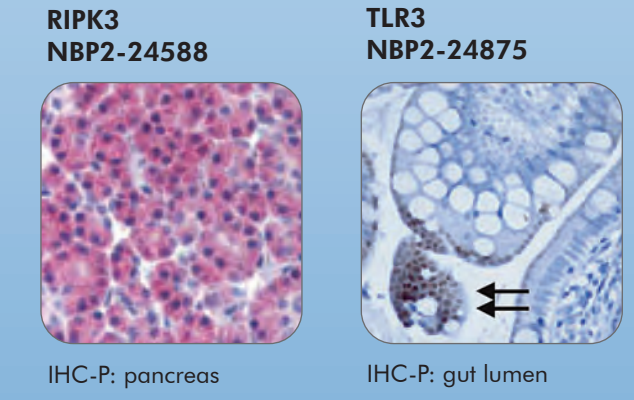
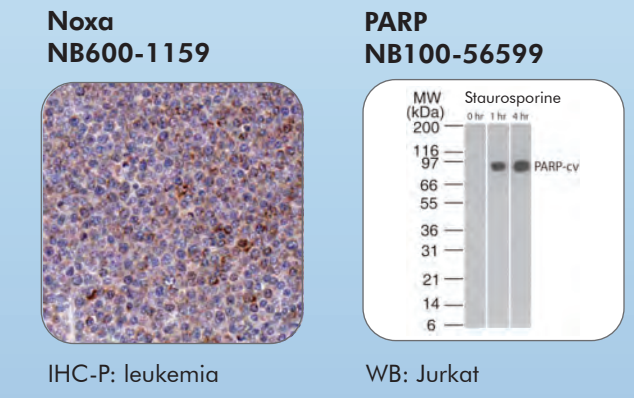
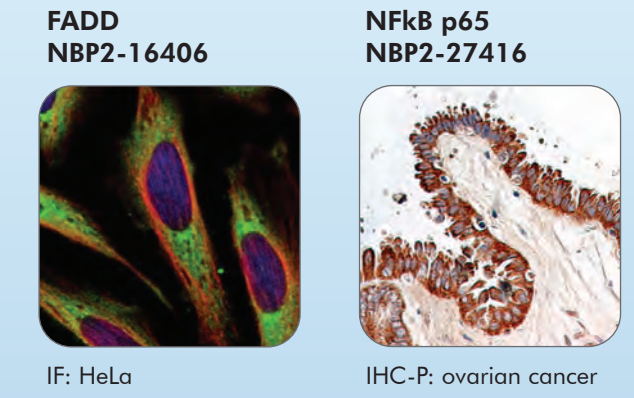
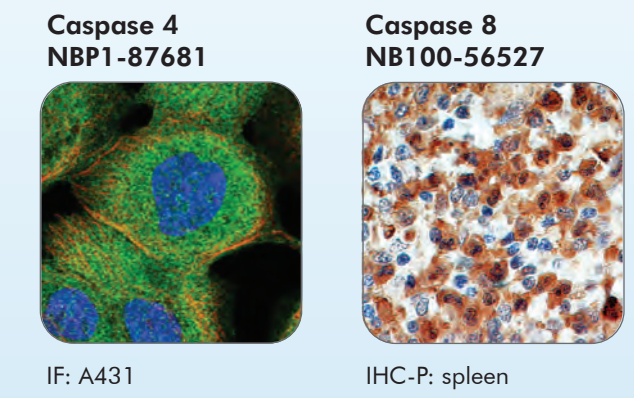
Autophagy

Autophagy is a basic catabolic mechanism involving lysosomal degradation and recycling of surplus or dysfunctional cellular components. Autophagy may promote cell survival or cell death depending on its execution context.



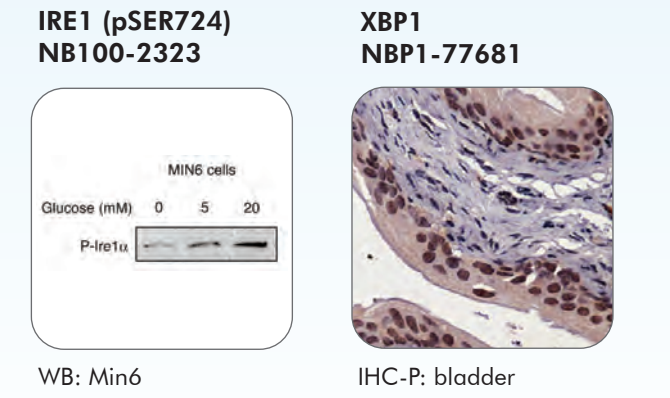
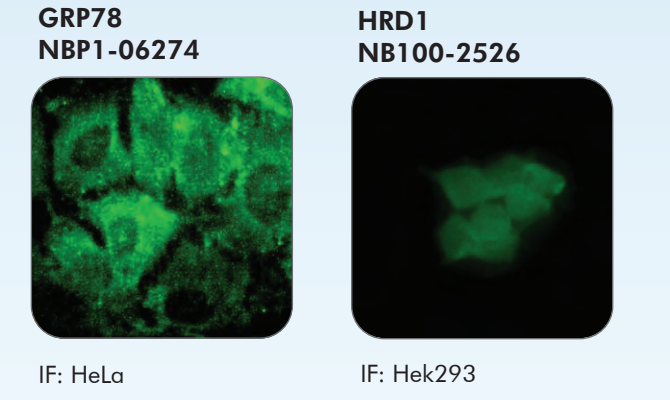
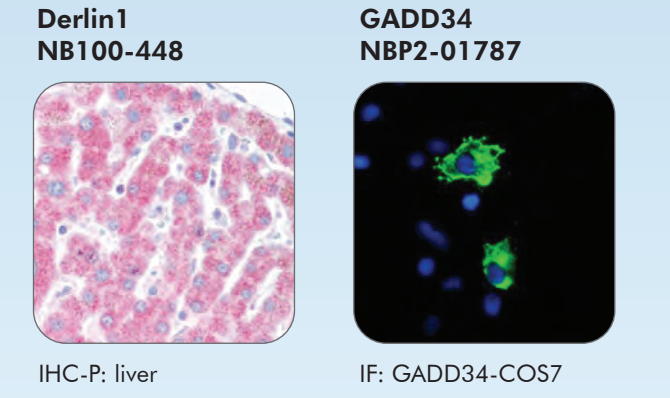
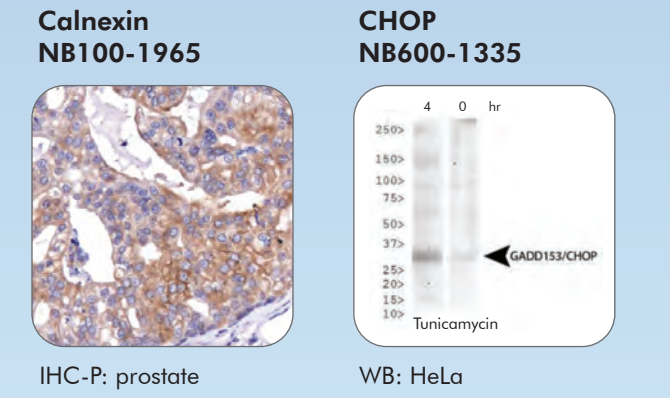
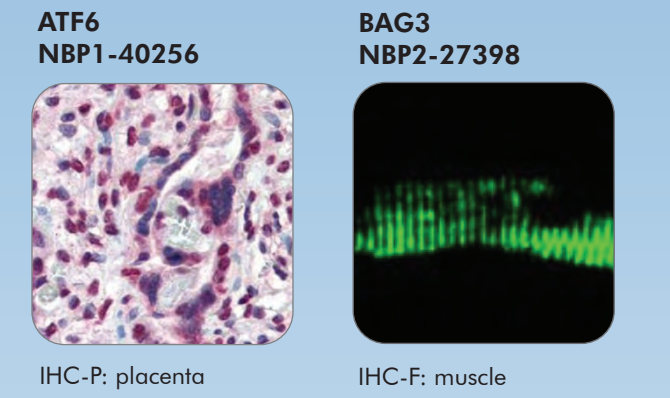
Necroptosis

Necroptosis is the execution of necrosis, a form of fatal cell injury, through regulated cell signaling pathways. Necroptosis plays a role in various pathologies, particularly those involving inflammatory processes.



UPR (Unfolded Protein Response)

The UPR is a stress response activated in response to an accumulation of unfolded or misfolded proteins in the endoplasmic reticulum. Active UPR initiates signaling pathways to either correct the problem or promote apoptosis.



Viability Dyes

7-AAD
Calcein Blue, AM
Calcein, AM
DAPI
Propidium Iodide

NB22-29446
NB22-31158
NB22-31157
NB22-31156
NB22-31155

Extraction Kits

Endoplasmic Reticulum Enrichment Kit
Mitochondrial Extraction Kit
Nuclear Extraction Kit

NB22-29482
NB22-29448
NB22-29447

Cell Death Analysis Kits

Annexin V-FITC Kit [FITC]
APO-BRDU (TUNEL) Kit
APO-BRDU-IHC (TUNEL) Kit
APO-DIRECT (TUNEL) Kit
Caspase Kit [FAM-VAD-FMK]
Caspase Kit [SR-VAD-FMK]
Mitochondrial Membrane Potential Kit
NF-kB p65 ELISA Kit
pSIVA™ Flow Kit [IANBD]
pSIVA™ Microscopy Kit [IANBD]

NB22-29373
NB22-31161
NB22-31164
NB22-31159
NB22-29383
NB22-29399
NB22-31093
NB22-31042
NB22-29611
NB22-29382

U.S. AND INTERNATIONAL CUSTOMERS

Phone: 303.730.1950
888.506.6887
Fax: +44 (0)871 971 1635
Email: orders@novusbio.com
Web: www.novusbio.com

CANADIAN CUSTOMERS

Phone: 905.827.6400
855.668.8722
Fax: 905.827.6402
Email: canada@novusbio.com
Web: www.novusbio.com/canada

EUROPEAN CUSTOMERS

NOVUS EUROPE
Phone: +44 (0)1223 426001
Fax: +44 (0)871 971 1635
Email: europe@novusbio.com

NOVUS BELGIUM
Phone: +32 2 401 22 53
Fax: +32 2 401 22 54
Email: belgium@novusbio.com

NOVUS DENMARK
Phone: +45 36 92 78 17
Fax: +45 36 92 78 18
Email: denmark@novusbio.com

NOVUS FINLAND
Phone: +358 9 231 954 48
Fax: +358 9 231 954 49
Email: finland@novusbio.com

NOVUS FRANCE
Phone: +33 1 76 77 45 30
Fax: +33 1 76 77 45 31
Email: france@novusbio.com

NOVUS GERMANY
Phone: +49 0800 723 5208
Fax: +49 0800 58926 79
Email: germany@novusbio.com

NOVUS ITALY
Phone: +39 02 4032 6786
Fax: +39 02 4032 6340
Email: italy@novusbio.com

NOVUS IRELAND
Phone: +353 1 506 0361
Fax: +353 1 506 0362
Email: ireland@novusbio.com

NOVUS NETHERLANDS
Phone: +31 2 07168336
Fax: +31 2 07168337
Email: netherlands@novusbio.com

NOVUS NORWAY
Phone: +47 21 03 42 86
Fax: +47 21 03 42 87
Email: norway@novusbio.com

NOVUS SWEDEN
Phone: +46 856619332
Fax: +46 856619333
Email: sweden@novusbio.com