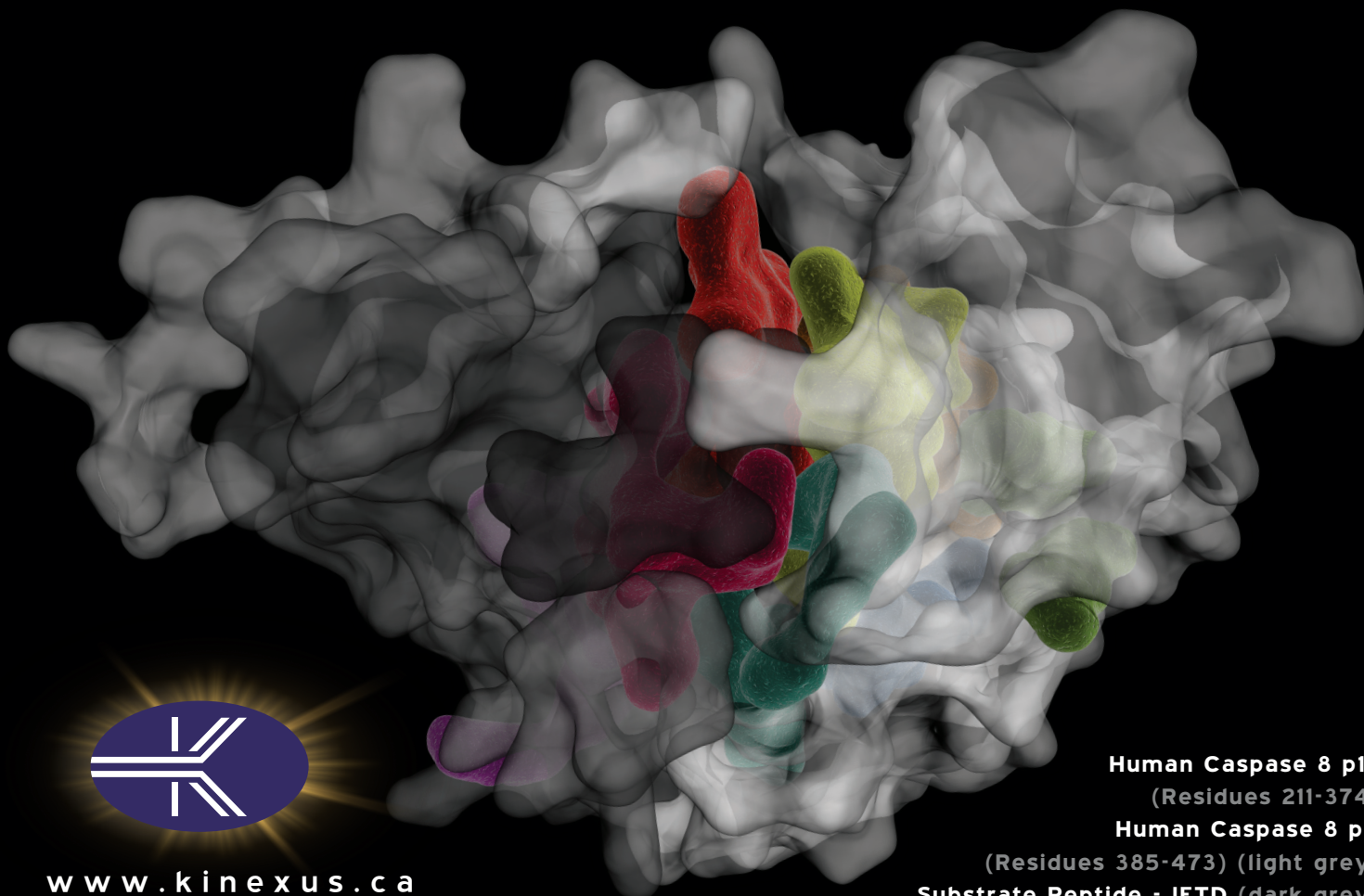




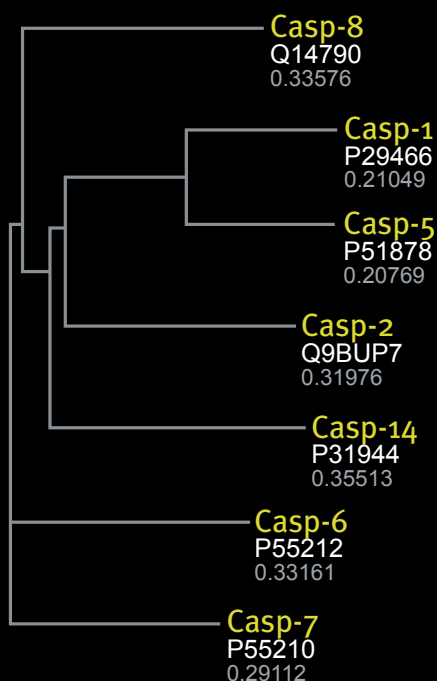
www.kinexus.ca

1-866-kinexus TOLL FREE PHONE

Human Caspase 8 p18
(Residues 211-374)

Human Caspase 8 p11
(Residues 385-473) (light grey)
Substrate Peptide - IETD (dark grey)
PDB: 1QTN

Caspase Catalytic Domains



Casp-8	235	YCLIIINNHNFAKAREKVPKLHSIRDNRNGTHLDAGALTTTFEELHFEIKPHDDCTVEQIYEILKIY	
Casp-1	164	LALIICNEEFDS-----IPRRTGAEVDTITGMTMLLQNLGYSDVVKKNLTASDMTTELEAF	
Casp-2	201	LALVLSNVHFTGE-----KELEFRSGGDVDHSTLVTLFKLLGYDVHVLCDQTAQEMQEKLNQF	
Casp-14	19	LALILCVT-----KAREGSEEDLDALEHMFRLRFESTMKRDPTAEQFQEELEKF	
Casp-6	46	IALIFNHERFFWH-----LTLPERRGTCADRDNLTTRFSDLGFVEKCFNDLKAEELLLKIHEV	
Casp-7	69	KCIIINNKNFDKV-----TGMGVRNGTDKDAEALFKCFRSLGFDVIVYNDSCSAKMQDLLKKA	
Casp-8	300	QLMDHSN--MDFCICILSHGDKGIITYGTGQ----EAPIYELTSQFTGLKCPSLAGK	
Casp-1	219	AHRPEHK--TSDSTFLVFMESHGIREGICGKKHSEQVPDILQLNAIFNMLNTKNCPSLKDK	
Casp-2	259	AQLPAHR--VTDSCIVAALSHGVEGATYGVGDG----KLLQLQEVFQLFDNANCPSLQNK	
Casp-14	69	QQAIDSREDPVSCAFVVLMAHGREGFLKGEDGE----MVKLENLFEALNNKNCQALRAK	
Casp-6	104	STVSHAD---ADCFVCVFLSHGEGNHIYAYDA-----KIEIQTLTGFLKGDKCHSLVGK	
Casp-7	127	SEEDHTN---AACFACILLSHGEEENVYIGKDG-----VTPIKDLTAHFRGDRCKTLLEK	
Casp-8	352	PKVFFIQACQGDNYQKG-----IPVETDSEEQPYLEMDLSSPQTRYIPDEADFL	
Casp-1	277	PKVIIQACRGD--SPGVVWFKDSVGVSGNLSLPTTEEFEDDAIKK-AHIEKDFIA	
Casp-2	312	PKMFFIQACRGDETDRGVDQDQGNHAGSPGCEESDAGKEKLPKMRLPTRSDMIC	
Casp-14	124	PKVYIIQACRG-----EQRDPGETVGGDEIVMVIKDSPTIPTYTDALH	
Casp-6	155	PKIFIIQACRGNQHDVPV--IPLDVVDNQTEKLDTNITEVDAASVYTL PAGADFLM	
Casp-7	178	PKLFFIQACRG-----TELDGDIQADSGPINDTDANPRYKIPVEADFLF	
Casp-8	402	GMATVNNCVSYRNPAEGTWYIQSLCQSLRERCPRGGDILLTEV	440
Casp-1	330	FCSSTPDNVSWRHPTMGSVFIGRLIEHMQEYA-CSCDVEEIFRKV	373
Casp-2	367	GYACLKGTAAAMRNTKRGSWYIEALAQVFSERA-CDMHVADMLVKV	410
Casp-14	168	VYSTVEGYIAYRHDQKGSCFIQTLVDVFTKRK---GHILELLTEV	209
Casp-6	209	CYSAEGYYSHRETVNGSWYIQDLCEMLGKYG--SSLEFTELLTLV	252
Casp-7	222	AYSTVPGYYSWRSPGRGSWFVQALCSILEEHG-KDLEIMQILTRV	265