

Caspase Catalytic Domain Alignment

Casp-8	235	YCLIIINNHNFAKAREKVPKLSIRDNRNGTHLDAGALTTTFFELHFEIKPHDDCTVEQIYEILKIY
Casp-1	164	LALIICNEEFDS-----IPRRTGAEVDITGMTLLQNLGYSVDVKKNLTASDMTTELEAF
Casp-2	201	LALVLSNVHFTGE-----KELEFRSGGDVDHSTLVTLFKLLGYDVHVLCDQTAQEMQEKLQNF
Casp-14	19	LALILCVT-----KAREGSEEDLDAL EHMFRQLRFESTMKRDPTAEQFQEELEKF
Casp-6	46	IALIFNHERFFWH-----LTLPERRGTCADRDNLTRRFSDLGFVKKCFNDLKAPELLLKIHEV
Casp-7	69	KCIIINNKNFDKV-----TGMGVRNGTDKDAEALFKCFRSLGFDVIVYND CSCAKMQDLLKKA
Casp-8	300	QLMDHSN---MDCFICILSHGDKGIYGTGQ-----EAPIYELTSQFTGLKCPFLAGK
Casp-1	219	AHRPEHK--TSDSTFLVFMHSHGIREGICGKKHSEQVPDILQLNAIFNMLNTKNCPSLKD
Casp-2	259	AQLPAHR--VTDSCIVAALSHGVEGAIYGVVG-----KLLQLQEVFQLFDNANCPQLQNK
Casp-14	69	QQAIDSREDPVSCAFVVLMAHGREGFLKGEDGE-----MVKLENLFEALNNKNCQALRAK
Casp-6	104	STVSHAD---ADCFCVFLSHGEGNHIYAYDA-----KIEIQTLTGLFKGDKCHSLVVGK
Casp-7	127	SEEDHTN--AACFACILLSHGEEVIYGKDG-----VTPIKDLTAHFRGDRCKTLLEK
Casp-8	352	PKVFFIQACQGDNYQKG-----IPVETDSEEQPYLEMDLSSPQTRYIPDEADFL
Casp-1	277	PKVIIIQACRGD-SPGVVWFKDSVGVSGNLSLPTTEEFEDDAIKK-AHIEKDFIA
Casp-2	312	PKMFFIQACRGDETDGVDQDQGNHAGSPGCEESDAGKEKLPKMRLPTRSDMIC
Casp-14	124	PKVYIIQACRG-----EQRDPGETVGGDEIVMVIKDSPTIPTYTALH
Casp-6	155	PKIFIIQACRGNQHDPVP-IPLDVVDNQTEKLDNITEVDAASVYTL PAGADFLM
Casp-7	178	PKLFFIQACRG-----TELDDGIQADSGPINDTDANPRYKIPVEADFLF
Casp-8	402	GMATVNNCVSYRNPAEGTWYIQSLCQSLRERCPRGDDILTILTEV 440
Casp-1	330	FCSSTPDNVSWRHPTMGSVFIGRLIEHMQEYA-CSCDVEEIFRKV 373
Casp-2	367	GYACLKGTAAAMRNTKRGSWYIEALAQVFSERA-CDMHVADMLVKV 410
Casp-14	168	VYSTVEGYIAYRHDQKGSFCFIQTLVDVFTKRK---GHILELLTEV 209
Casp-6	209	CYSVAEGYYSHRETVNGSWYIQDLCEMLGKYG-SSLEFTELLTLV 252
Casp-7	222	AYSTVPGYYSWRSPGRGSWFVQALCSILEEHG-KDLEIMQILTRV 265

Human Caspase 8 p18
(residues 211-374, white)

Human Caspase 8 p11
(residues 385-473, light
grey)

Substrate Peptide IETD
(dark grey)

PDB: 1QTN

