

Dipeptide Library

The table below presents all dipeptides possibilities. Dipeptides containing N-terminal Glutamine or Glutamic acid are known to form pGlu. The purity of the grey highlighted peptides is expressed as a sum of both products (pGlu formation can vary from 20% to 50%), which is >95% pure by analytical HPLC. Once dissolved in solutions, these dipeptides should be utilized fairly quickly. Empty boxes denote that the peptides are not available. Unless otherwise specified, all other peptides are >95%.

Dipeptides

AA	DA	EA	FA	GA	HA	IA	KA	LA	MA	NA	PA	QA	RA	SA	TA	VA	WA	YA
AD	DD	ED	FD	GD	HD	ID	KD	LD	MD	ND	PD	QD	RD	SD	TD	VD	WD	YD
AE	-	EE	FE	GE	HE	IE	KE	LE	ME	NE	PE	QE	RE	SE	TE	VE	WE	YE
AF	DF	-	FF	GF	HF	IF	KF	LF	MF	NF	PF	QF	RF	SF	TF	VF	WF	YF
AG	DG	EG	FG	GG	-	-	KG	LG	MG	NG	PG	QG	RG	-	TG	VG	WG	YG
AH	DH	EH	FH	GH	HH	IH	KH	LH	MH	NH	PH	QH	RH	SH	TH	VH	WH	YH
AI	-	EI	FI	GI	HI	II	KI	LI	MI	NI	PI	QI	RI	SI	TI	VI	WI	YI
AK	DK	EK	FK	-	HK	IK	KK	-	MK	NK	PK	QK	RK	SK	TK	VK	WK	YK
AL	DL	EL	FL	-	HL	IL	KL	LL	ML	NL	-	QL	RL	SL	TL	VL	WL	YL
AM	DM	EM	FM	GM	HM	IM	KM	LM	MM	NM	PM	QM	RM	SM	TM	VM	WM	YM
-	DN	EN	FN	GN	HN	IN	KN	LN	MN	NN	PN	QN	RN	SN	TN	VN	WN	YN
AP	DP	EP	-	GP	HP	IP	KP	-	-	NP	PP	QP	RP	SP	TP	VP	WP	YP
AQ	DQ	EQ	FQ	GQ	-	IQ	KQ	-	MQ	NQ	PQ	QQ	-	SQ	TQ	VQ	WQ	YQ
AR	DR	ER	FR	GR	HR	IR	KR	-	MR	NR	-	-	RR	SR	TR	VR	WR	YR
AS	DS	ES	FS	GS	HS	IS	KS	-	MS	-	PS	QS	RS	SS	TS	VS	WS	YS
AT	DT	ET	FT	GT	HT	IT	KT	LT	MT	NT	PT	QT	RT	ST	TT	VT	WT	YT
AV	DV	EV	FV	GV	HV	IV	KV	LV	MV	NV	PV	QV	RV	SV	TV	-	VV	YV
AW	DW	EW	FW	GW	HW	IW	KW	LW	MW	NW	PW	QW	RW	SW	TW	VW	WW	YW
AY	DY	EY	FY	GY	HY	-	KY	LY	MY	NY	PY	QY	-	SY	TY	VY	WY	YY