

IDseek® STR Panel — Loci Reference: CODIS & ENFSI Designations

OmniSTR Global • mYSTR • CombiSTR Plus — NimaGen, distributed by Promega Corporation (US)

Locus	Designation	OmniSTR	mYSTR	CombiSTR+		Locus	Designation	OmniSTR	mYSTR	CombiSTR+
AUTOSOMAL STR LOCI						Y-CHROMOSOMAL STR LOCI				
Amelogenin	Sex marker	✓	—	✓		DYS19	^ ENFSI	✓	✓	✓
CSF1PO	# CODIS+ENFSI	✓	—	✓		DYS385a/b	^ ENFSI	—	✓	✓
D1S1656	# CODIS+ENFSI	✓	—	✓		DYF387S1a/b	Additional	—	✓	✓
D2S441	# CODIS+ENFSI	✓	—	✓		DYS389(I/II)	^ ENFSI	—	✓	✓
D2S1338	# CODIS+ENFSI	✓	—	✓		DYS390	^ ENFSI	—	✓	✓
D3S1358	# CODIS+ENFSI	✓	—	✓		DYS391	^ ENFSI	✓	✓	✓
D4S2408	^ ENFSI	✓	—	✓		DYS392	^ ENFSI	—	✓	✓
D5S818	# CODIS+ENFSI	✓	—	✓		DYS393	^ ENFSI	—	✓	✓
D6S1043	^ ENFSI	✓	—	✓		DYS437	^ ENFSI	—	✓	✓
D7S820	# CODIS+ENFSI	✓	—	✓		DYS438	^ ENFSI	—	✓	✓
D8S1179	# CODIS+ENFSI	✓	—	✓		DYS439	^ ENFSI	—	✓	✓
D9S1122	^ ENFSI	✓	—	✓		DYS448	^ ENFSI	—	✓	✓
D10S1248	# CODIS+ENFSI	✓	—	✓		DYS449	Additional	—	✓	✓
D12S391	# CODIS+ENFSI	✓	—	✓		DYS456	^ ENFSI	—	✓	✓
D13S317	# CODIS+ENFSI	✓	—	✓		DYS458	^ ENFSI	—	✓	✓
D16S539	# CODIS+ENFSI	✓	—	✓		DYS460	Additional	—	✓	✓
D17S1301	^ ENFSI	✓	—	✓		DYS461	Additional	—	✓	✓
D18S51	# CODIS+ENFSI	✓	—	✓		DYS481	Additional	—	✓	✓
D19S433	# CODIS+ENFSI	✓	—	✓		DYS505	Additional	—	✓	✓
D20S482	^ ENFSI	✓	—	✓		DYS518	Additional	—	✓	✓
D21S11	# CODIS+ENFSI	✓	—	✓		DYS522	Additional	—	✓	✓
D22S1045	# CODIS+ENFSI	✓	—	✓		DYS533	Additional	—	✓	✓
FGA	# CODIS+ENFSI	✓	—	✓		DYS549	Additional	—	✓	✓
Penta D	Additional	✓	—	✓		DYS570	Additional	—	✓	✓
Penta E	Additional	✓	—	✓		DYS576	Additional	—	✓	✓
SE33	^ ENFSI	✓	—	✓		DYS612	Additional	—	✓	✓
TH01	# CODIS+ENFSI	✓	—	✓		DYS627	Additional	—	✓	✓
TPOX	# CODIS+ENFSI	✓	—	✓		DYS635	^ ENFSI	—	✓	✓
vWA	# CODIS+ENFSI	✓	—	✓		DYS643	Additional	—	✓	✓
						Y-GATA-H4	^ ENFSI	—	✓	✓
						TXLNKY	Sex marker	—	✓	✓

LEGEND:

CODIS + ENFSI

^ ENFSI only

Additional

Sex marker

✓ = Present in kit

— = Absent

OmniSTR: 29 mYSTR: 31 CombiSTR Plus: 60 loci