

# Finnzymes to Bio-Rad Reagent Conversion Chart

MJR Catalog No.	MJR Product Description	Equivalent Bio-Rad Product	Bio-Rad Catalog No.
F-410S F-410L	DyNAmo hot start SYBR Green qPCR kit, 40 rxns DyNAmo hot start SYBR Green qPCR kit, 200 rxns		
F-400S F-400L	DyNAmo SYBR Green qPCR kit, 40 rxns DyNAmo SYBR Green qPCR kit, 200 rxns	iQ SYBR Green supermix or iTaq SYBR Green supermix with ROX	170-8880, 170-8882, 170-8888, 170-8885 or 170-8850, 170-8851, 170-8852, 170-8853
F-400RS F-400RL	DyNAmo SYBR Green qPCR kit with ROX, 40 rxns DyNAmo SYBR Green qPCR kit with ROX, 200 rxns		
F-420S F-420L	DyNAmo Capillary kit, 100 rxns DyNAmo Capillary kit, 500 rxns	No equivalent Bio-Rad product; please contact Finnzymes at <a href="http://www.finnzymes.fi">www.finnzymes.fi</a>	
F-500S F-500L F-502S F-502L	Native DyNAzyme DNA polymerase, 100 U Native DyNAzyme DNA polymerase, 500 U Native DyNAzyme DNA polymerase (Mg-free buffer), 100 U Native DyNAzyme DNA polymerase (Mg-free buffer), 500 U	iTaq hot-start DNA polymerase	170-8870 or 170-8875
F-501L F-501S F-503L F-503S	Recombinant DyNAzyme II DNA polymerase, 250 U Recombinant DyNAzyme II DNA polymerase, 1,000 U Recombinant DyNAzyme II DNA polymerase (Mg-free buffer), 250 U Recombinant DyNAzyme II DNA polymerase (Mg-free buffer), 1,000 U		
F-505S F-505L	DyNAzyme EXT, 200 U DyNAzyme EXT, 1,000 U	iProof high-fidelity DNA polymerase	172-5300, 172-5301 and 172-5302
F-550S F-550L F-551S F-551L F-552XS F-552S F-552L	Native DyNAzyme DNA polymerase kit, 100 U Native DyNAzyme DNA polymerase kit, 500 U Recombinant DyNAzyme II DNA polymerase kit, 100 U Recombinant DyNAzyme II DNA polymerase kit, 500 U DyNAzyme EXT DNA polymerase kit, 50 U DyNAzyme EXT DNA polymerase kit, 200 U DyNAzyme EXT DNA polymerase kit, 1000 U	iTaq hot-start DNA polymerase and dNTP mix (please note: Bio-Rad does not yet offer iTaq DNA polymerase in a kit along with control templates, primers, and mass standards)	170-8870 or 170-8875 and 170-8874
F-530 F-530S F-530L	Phusion high-fidelity DNA polymerase, 20 U Phusion high-fidelity DNA polymerase, 100 U Phusion high-fidelity DNA polymerase, 500 U	iProof high-fidelity DNA polymerase	172-5300, 172-5301 and 172-5302
F-531S F-531L	Phusion master mix with high-fidelity buffer, 100 rxns Phusion master mix with high-fidelity buffer, 500 rxns	iProof master mix with high-fidelity buffer	172-5310 and 172-5311
F-532S F-532L	Phusion master mix with GC buffer, 100 rxns Phusion master mix with GC buffer, 500 rxns	iProof master mix with GC buffer	172-5320 and 172-5321
F-553S F-553L	Phusion high-fidelity DNA polymerase kit, 50 U Phusion high-fidelity DNA polymerase kit, 200 U	iProof high-fidelity DNA polymerase kit	172-5330 and 172-5331
F-560S F-560L F-560XL	dNTP mix solution, 100 rxns dNTP mix solution, 500 rxns dNTP mix solution, 2,500 rxns	dNTP mix	170-8874
F-570S F-570L F-580S F-580L	AMV reverse transcriptase, 500 U AMV reverse transcriptase, 2,500 U RobusT RT-PCR kit, 20 rxns RobusT RT-PCR kit, 100 rxns	iScript cDNA synthesis kit iScript Select cDNA synthesis kit or iScript one-step quantitative RT-PCR kits	170-8890 and 170-8891, or 170-8896 and 170-8897, or 170-8892, 170-8893, 170-8894 and 170-8895



# Finnzymes to Bio-Rad Reagent Conversion Chart

MJR Catalog No.	MJR Product Description	Equivalent Bio-Rad Product	Bio-Rad Catalog No.
F-700	Template generation system kit, 20 rxns		
F-702	Template generation system II kit, 20 rxns		
F-778	Entranceposon (camR), 20 rxns	No equivalent Bio-Rad product; please contact Finnzymes at <a href="http://www.finnzymes.fi">www.finnzymes.fi</a>	
F-779	Entranceposon (kanR), 20 rxns		
F-784	Entranceposon (tetR), 20 rxns		
F-701	Mutation generation system kit, 10 rxns		
F-750	MuA transposase with 5x buffer, 20 rxns	No equivalent Bio-Rad product; please contact Finnzymes at <a href="http://www.finnzymes.fi">www.finnzymes.fi</a>	
F-750C	MuA transposase (5x concentrate), 20 rxns		
F-760	Entranceposon (M1-camR) 20 rxns		
F-762	Entranceposon (M1-kanR) 20 rxns		
F-703	STOP Generation Kit, 10 rxns	No equivalent Bio-Rad product; please contact Finnzymes at <a href="http://www.finnzymes.fi">www.finnzymes.fi</a>	
F-751	Entranceposon, (camR) 20 rxns	No equivalent Bio-Rad product; please contact Finnzymes at <a href="http://www.finnzymes.fi">www.finnzymes.fi</a>	
F-759	Entranceposon, (kanR) 20 rxns		
F-764	Entranceposon, (tetR) 20 rxns		
F-765	pEntranceposon, (camR) plasmid		
F-766	pEntranceposon, (kanR) plasmid		
F-767	pEntranceposon, (tetR) plasmid		
F-768	Entranceposon, (camR-2) 20 rxns		
F-771	Entranceposon, (SupF) 20 rxns		
F-773	pEntranceposon, (SupF) plasmid		
F-774	Entranceposon, (lacZ) 20 rxns		
F-775	Entranceposon, (lacZ PT) 20 rxns		
F-800S	Ampli-Y bovine embryo sexing kit, 25 rxns	No equivalent Bio-Rad product; please contact Finnzymes at <a href="http://www.finnzymes.fi">www.finnzymes.fi</a>	
F-800L	Ampli-Y bovine embryo sexing kit, 100 rxns		
MJR Item No.	Accessory Items	Equivalent Bio-Rad Product	
F-301L	Lambda-HindIII, 500 ug		
F-302LD	φX-HaeIII, 100 ug with dye		
F-302SD	φX-HaeIII, 20 ug with dye		
F-303LD	DNA marker mix, large		
F-303SD	DNA marker mix, small		
F-324-1	DyNAfreezer, 24 X 1.5 ml		
F-332-1	DyNAfreezer, 12 X 1.5 ml		
F-340	CapEase tool, blue base and white lid		
F-341	CapEase tool, transparent		
F-510L	Mg-free DyNAzyme buffer with Mg solution		
F-511L	Optimized DyNAzyme buffer		
F-512L	Mg-free DyNAzyme EXT buffer with Mg solution		
F-514L	Optimized DyNAzyme EXT buffer	No equivalent Bio-Rad product; please contact Finnzymes at <a href="http://www.finnzymes.fi">www.finnzymes.fi</a>	
F-516L	DyNAzyme detergent-free buffer		
F-517L	DyNAzyme EXT detergent-free buffer		
F-518L	Phusion high-fidelity buffer, 3 ml		
F-518S	Phusion high-fidelity buffer, 6 ml		
F-519	Phusion GC buffer, 1.5 ml		
F-519L	Phusion GC buffer, 6 ml		
F-519S	Phusion GC buffer, 3 ml		
F-520L	HPLC Phusion high-fidelity buffer, 3 ml		
F-520S	HPLC Phusion high-fidelity buffer, 6 ml		
F-521L	HPLC Phusion GC buffer, 6 ml		
F-521S	HPLC Phusion GC buffer, 3 ml		
F-600	DyNAwax, 20 ml		