

Flexi® Vector Name <sup>1</sup>	Cat. #	Drug Selection <sup>2</sup>	Expression Application	Promoter for Expression			Peptide Fusion Tag	
				<i>E. coli</i>	Mammalian Cells	Cell-Free Transcription/ Translation	N-terminus	C-terminus
pF1A T7 Flexi® Vector pF1K T7 Flexi® Vector	C8441 C8451	Ampicillin Kanamycin	Inducible expression of native protein	T7		T7		
pFN2A (GST) Flexi® Vector pFN2K (GST) Flexi® Vector	C8461 C8471	Ampicillin Kanamycin	Soluble expression and purification via an N-terminal GST tag	T7		T7	GST	
pF3A WG (BYDV) Flexi® Vector pF3K WG (BYDV) Flexi® Vector	L5671 L5681	Ampicillin Kanamycin	In vitro wheat germ expression of native protein			T7, SP6		
pF25A ICE T7 Flexi® Vector pF25K ICE T7 Flexi® Vector	L1061 L1081	Ampicillin Kanamycin	In vitro insect cell free expression of native protein			T7		
pFN6A (HQ) Flexi® Vector pFN6K (HQ) Flexi® Vector	C8511 C8521	Ampicillin Kanamycin	Inducible expression and protein purification via a metal affinity resin	T7		T7	HQHQHQ	
pFC7A (HQ) Flexi® Vector pFC7K (HQ) Flexi® Vector	C8531 C8541	Ampicillin Kanamycin	Inducible expression and protein purification via a metal affinity resin	T7		T7		HQHQHQ
pF4A CMV Flexi® Vector pF4K CMV Flexi® Vector	C8481 C8491	Ampicillin Kanamycin	Constitutive high expression of native protein		CMV	T7		
pF5A CMV-neo Flexi® Vector pF5K CMV-neo Flexi® Vector	C9401 C9411	Ampicillin/Neomycin Kanamycin/Neomycin <sup>M</sup>	Constitutive high expression of native protein with selection for stable transfectants		CMV	T7		
pF9A CMV <i>hRluc</i> -neo Flexi® Vector	C9361	Ampicillin Neomycin <sup>M</sup>	Constitutive high expression of native protein with selection and reporter screening for stable transfectants		CMV	T7		
pFN10A (ACT) Flexi® Vector	C9331	Ampicillin/Neomycin	Mammalian in vivo protein protein interaction		CMV	T7	HSV VP16 activation domain	
pN11A (BIND) Flexi® Vector	C9341	Ampicillin	Mammalian in vivo protein: protein interaction		CMV	T7	GAL4 DNA-binding domain	
pF12A RM Flexi® Vector pF12K RM Flexi® Vector	C9431 C9441	Ampicillin Kanamycin	Regulated mammalian protein expression		12λOP-Mini CMV			

<sup>1</sup> The "pF" indicates that the vector is a Flexi® Vector. The letter after "pF" indicates the position of any expression tags ("N" for an N-terminal expression tag and "C" for a C-terminal expression tag). The number associated with the vector specifies the type of expression and application. The letters "A" and "K" designate the bacterial drug selection for the vector ("A"=ampicillin and "K"=kanamycin).

<sup>2</sup> The "M" indicates the vector provides resistance to the indicated drug in mammalian cells.

Flexi® Vector Name <sup>1</sup>	Cat.#	Drug Selection <sup>2</sup>	Expression Application	Promoter for Expression			Peptide Fusion Tag	
				<i>E. coli</i>	Mammalian Cells	Cell-Free Transcription/ Translation	N-terminus	C-terminus
pFN18A HaloTag® T7 Flexi® Vector pFN18K HaloTag® T7 Flexi® Vector	G2751 G2681	Ampicillin Kanamycin	Protein interaction analysis and protein purification	T7		T7	HaloTag®	
pFN19A HaloTag® T7 SP6 Flexi® Vector pFN19K HaloTag® T7 SP6 Flexi® Vector	G1891 G1841	Ampicillin Kanamycin	Protein interaction analysis	T7		T7, SP6	HaloTag®	
pFC20A HaloTag® T7 SP6 Flexi® Vector pFC20K HaloTag® T7 SP6 Flexi® Vector	G1681 G1691	Ampicillin Kanamycin	Protein interaction analysis	T7		T7, SP6		HaloTag®
pFC14A HaloTag® CMV Flexi® Vector pFC14K HaloTag® CMV Flexi® Vector	G9651 G9661	Ampicillin Kanamycin	Cell imaging and protein interaction analysis; high constitutive expression		CMV			HaloTag®
pFC15A HaloTag® CMVd1 Flexi® Vector pFC15K HaloTag® CMVd1 Flexi® Vector	G1611 G1601	Ampicillin Kanamycin	Cell imaging and protein interaction analysis; medium constitutive expression	T7	CMVd1	T7, SP6		HaloTag®
pFC16A HaloTag® CMVd2 Flexi® Vector pFC16K HaloTag® CMVd2 Flexi® Vector	G1591 G1571	Ampicillin Kanamycin	Cell imaging and protein interaction analysis; low constitutive expression	T7	CMVd2	T7, SP6		HaloTag®
pFC17A HaloTag® CMVd3 Flexi® Vector pFC17K HaloTag® CMVd3 Flexi® Vector	G1551 G1321	Ampicillin Kanamycin	Cell imaging and protein interaction analysis; very low constitutive expression	T7	CMVd3	T7, SP6		HaloTag®
pFN21A HaloTag® CMV Flexi® Vector pFN21K HaloTag® CMV Flexi® Vector	G2821 G2831	Ampicillin Kanamycin	Cell imaging and protein interaction analysis; high constitutive expression		CMV	T7	HaloTag®	
pFN22A HaloTag® CMVd1 Flexi® Vector pFN22K HaloTag® CMVd1 Flexi® Vector	G2841 G2851	Ampicillin Kanamycin	Cell imaging and protein interaction analysis; medium constitutive expression	T7	CMVd1	T7, SP6	HaloTag®	
pFN23A HaloTag® CMVd2 Flexi® Vector pFN23K HaloTag® CMVd2 Flexi® Vector	G2861 G2871	Ampicillin Kanamycin	Cell imaging and protein interaction analysis; low constitutive expression	T7	CMVd2	T7, SP6	HaloTag®	
pFN24A HaloTag® CMVd3 Flexi® Vector pFN24K HaloTag® CMVd3 Flexi® Vector	G2881 G2981	Ampicillin Kanamycin	Cell imaging and protein interaction analysis; very low constitutive expression	T7	CMVd3	T7, SP6	HaloTag®	

<sup>1</sup> The "pF" indicates that the vector is a Flexi® Vector. The letter after "pF" indicates the position of any expression tags ("N" for an N-terminal expression tag and "C" for a C-terminal expression tag). The number associated with the vector specifies the type of expression and application. The letters "A" and "K" designate the bacterial drug selection for the vector ("A"=ampicillin and "K"=kanamycin).

<sup>2</sup> The "M" indicates the vector provides resistance to the indicated drug in mammalian cells.



AF9GE096 0706GE096



Promega Corporation • 2800 Woods Hollow Road • Madison, WI 53711-5399 USA • Telephone 608-274-4330 • Fax 608-277-2601  
www.promega.com

©2008 Promega Corporation. All Rights Reserved.

Prices and specifications subject to change without prior notice.

Printed in USA Rev. 05/09

17687-FL-GN

Part #GE096

