

Technical Bulletin

pAdVAntage[™] Vector

INSTRUCTIONS FOR USE OF PRODUCT E1711.

PRINTED IN USA. Revised 11/06

Part# TB207

vww.promega.com

pAdVAntage[™] Vector

All technical literature is available on the Internet at: www.promega.com/tbs/ Please visit the web site to verify that you are using the most current version of this Technical Bulletin. Please contact Promega Technical Services if you have questions on use of this system. E-mail: techserv@promega.com

I.	Description	1
II.	Product Components and Storage Conditions	2
III.	 General Considerations A. Optimization of the Ratio of pAdVAntage[™] Vector to Expression Vector DNA B. Sample Experiments Demonstrating pAdVAntage[™] Vector Enhancement of Expression in HeLa and 293 Cells 	4
IV.	pAdVAntage™ Vector Restriction Sites	5
v.	Related Products	7
VI.	References	7

I. Description

Co-transfection of mammalian cells with the pAdVAntage[™] Vector enhances transient protein expression in a variety of cell types by increasing translation initiation.

Transfection of mammalian cells with an expression vector often results in suboptimal expression of the protein of interest. Double-stranded RNA (dsRNA) generated during transfection is thought to activate the dsRNA-activated inhibitor (DAI), one of several enzymes involved in the host cell's antiviral defense system. DAI phosphorylates the translation initiation factor eIF-2, halting translation and therefore protein production (1, reviewed in reference 2; see Figure 1).

However, inhibition of translation by DAI can be overcome with the adenoviral Virus Associated I RNA (VAI RNA) that is produced by RNA polymerase III following co-transfection with the pAdVAntage[™] Vector. The VAI RNA binds to DAI, preventing its activation, thereby allowing translation and protein expression (3,4, reviewed in reference 2; see Figure 1).

The pAdVAntage[™] Vector contains base pairs 9,831–11,555 of the adenovirus type 2 genome on a 1,724bp SalI–HindIII fragment that encodes the virus-associated RNA genes, VAI and VAII.

 Promega Corporation
 2800
 Woods
 Hollow Road
 Madison, WI
 53711-5399
 USA

 Toll Free in USA 800-356-9526
 Phone 608-274-4330
 Fax
 608-277-2516
 www.promega.com

 Printed in USA.
 Part# TB207
 Part# TB207
 Page 1



II. Product Components and Storage Conditions

Product	Size	Cat.#
pAdVAntage™ Vector	20µg	E1711
pAdVAntage [™] Vector is supplied frozen in TE buffer.		

Storage Conditions: Store at -20°C.



Figure 1. Translation initiation and the effects of VAI RNA on this process.





Figure 2. pAdVAntage[™] Vector circle map and sequence reference points.

Sequence reference points:

Adenovirus DNA insert	1-1725
Adenovirus VAI RNA gene	780-939
Adenovirus VAII RNA gene	1036-1198
β-lactamase (Amp ^r) coding region	2907-3764

III. General Considerations

The level of expression enhancement attained by co-transfection with the pAdVAntageTM Vector is dependent upon a number of parameters, including the cell type, method of transfection and the ratio of pAdVAntageTM Vector DNA to expression vector DNA. Therefore, optimization of these parameters is critical for best results.

III.A. Optimization of the Ratio of pAdVAntage™ Vector to Expression Vector DNA

To optimally enhance protein expression following co-transfection with the pAdVAntage[™] Vector, experimentally determine the optimal ratio of pAdVAntage[™] Vector to expression vector DNA for each transfection system. For initial optimization experiments, we recommend using ratios of pAdVAntage[™] Vector to expression vector DNA in the range of 1:10 to 1:1, as demonstrated in the sample in Table 1.

	I	ONA Ratios	5	
(pAdVAn	tage™ Vecto	or DNA:Exp	pression Vec	tor DNA)
Transfected DNA	1:10	1:5	1:2.5	1:1
pAdVAntage™ Vector DNA (or pBR322 DNA in control plates)	0.5µg	1µg	2μg	5µg
Expression Vector DNA	5µg	5µg	5µg	5µg

Table 1. Sample Optimization of the Ratio of Co-Transfected DNAs.

III.B. Sample Experiments Demonstrating pAdVAntage™ Vector Enhancement of Expression in HeLa and 293 Cells

To determine how well the pAdVAntage[™] Vector can enhance protein expression, we performed co-transfection experiments of 293 and HeLa cells with pGL2-Control DNA and either pAdVAntage[™] DNA or pBR322 DNA as a control.

The cells were transfected in 60mm tissue culture dishes with 5µg of the pGL2-Control DNA (Cat.# E1611) using the ProFection® Mammalian Transfection System – Calcium Phosphate (Cat.# E1200). The ratio of pAdVAntage™ DNA or pBR322 DNA co-transfected with the pGL2-Control DNA ranged from 1:10 to 1:1 as detailed in Table 1. The cells were lysed forty-eight hours post-transfection, and luciferase activity expressed from the pGL2-Control DNA was measured using the Luciferase Assay System (Cat.# E1500). The relative-fold increase in luciferase activity due to co-transfection with the pAdVAntage™ Vector was determined using the luciferase activity of the control cells co-transfected with pBR322 as a baseline for comparison.

In these experiments, 293 cells co-transfected with the pAdVAntage[™] Vector demonstrated a 10- to 70-fold increase in luciferase expression from pGL2-Control DNA, while HeLa cells demonstrated a 4- to 40-fold increase in luciferase expression when co-transfected with the pAdVAntage[™] Vector.

IV. pAdVAntage[™] Vector Restriction Sites

The following restriction enzyme tables were constructed using DNASTAR[®] sequence analysis software. Please note that we have not verified this information by restriction digestion with each enzyme listed. The location given specifies the 3' end of the cut DNA (the base to the left of the cut site). For more information on the cut sites of these enzymes or if you identify a discrepancy, please contact your local Promega Branch or Distributor. In the U.S., contact Promega Technical Services at 800-356-9526. Vector sequences are also available from GenBank[®] database (GenBank[®]/EMBL Accession Number **U47294**) and on the Internet at: **www.promega.com/vectors/**

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AatII	2	914, 3899	CspI	1	1631
AccI	2	2, 314	Csp45I	1	841
AccIII	1	1171	DraI	3	2843, 2862, 3554
Acc65I	1	4372	DraII	3	1084, 1407, 3953
AflIII	2	1656, 2084	DrdI	2	2192, 4061
Alw44I	3	2398, 3644, 4141	DsaI	5	701, 790, 1325,
AlwNI	3	494, 1500, 2500			1363, 1489
ApaI	1	1377	EagI	3	855, 1108, 1664
AvaI	3	1101, 1528, 4376	EarI	5	215, 791, 1968,
BalI	2	382, 982			3772, 4260
BamHI	2	850, 4381	EclHKI	2	321, 2977
BbeI	3	207, 1371, 4203	Eco47III	1	930
BbrPI	1	1659	Eco52I	3	855, 1108, 1664
BbsI	1	566	Eco72I	1	1659
BglI	4	265, 726, 3097,	EcoICRI	2	425, 4368
-		4215	EcoNI	1	306
BsaI	1	3038	EcoRI	1	4360
BsaAI	2	1485, 1659	EcoRV	2	499, 595
BsaMI	1	44	EheI	3	205, 1369, 4201
BsmI	1	44	FseI	1	1110
Bsp120I	1	1373	FspI	5	43, 682, 734, 3199
BspHI	3	2804, 3812, 3917			4222
BspMI	2	77, 310	HincII	3	3, 398, 743
BssHII	4	646, 992, 1648,	HindII	3	3, 398, 743
		1650	HindIII	1	1725
BssSI	3	2257, 3641, 3948	HpaI	1	398
BstEII	3	407, 878, 1677	KasI	3	203, 1367, 4199
BstXI	1	21	KpnI	1	4376
BstZI	3	855, 1108, 1664	NaeI	2	629, 1108
Cfr10I	5	121, 627, 720,	NarI	3	204, 1368, 4200
		1106, 3057	NcoI	1	701

Table 1. Restriction Enzymes That Cut the pAdVAntage $^{\rm TM}$ Vector Between 1 and 5 Times.

Note: The enzymes listed in boldface type are available from Promega.

 Promega Corporation
 2800
 Woods
 Hollow Road
 Madison, WI
 53711-5399
 USA

 Toll Free in USA 800-356-9526
 Phone 608-274-4330
 Fax 608-277-2516
 www.promega.com

 Printed in USA.
 Part# TB207

 Revised 11/06
 Page 5

IV. pAdVAntage[™] Vector Restriction Sites (continued)

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
NdeI	1	4148	SacI	2	427, 4370
NgoMIV	2	627, 1106	SacII	3	1328, 1366, 1492
NheI	1	969	Sall	1	1
NotI	1	1664	ScaI	1	3457
NruI	1	1508	SmaI	1	4378
NsiI	1	1214	SplI	2	1481, 1685
NspI	3	1280, 2088, 4005	SspI	1	3781
PmlI	1	1659	StyI	4	20, 701, 1067,
Ppu10I	1	1210			1447
PpuMI	1	1084	TfiI	2	1919, 2059
Psp5II	1	1084	Tth111I	1	468
PspAI	1	4376	VspI	3	1855, 1914, 3149
PvuI	2	3347, 4243	XbaI	2	749, 4387
PvuII	3	1458, 1908, 4272	XmaI	1	4376
RsrII	1	1631	XmnI	1	3576

Table 1. Restriction Enzymes That Cut the pAdVAntage $^{\rm TM}$ Vector Between 1 and 5 Times (continued).

	Table 2.	Restriction	Enzymes	That Do	Not Cut the	pAdVAntage™	Vector.
--	----------	-------------	---------	---------	-------------	-------------	---------

AccB7I	BlpI	DraIII	PmeI	SphI
AfIII	Bpu1102I	Eco81I	PshAI	SrfI
AgeI	BsaBI	I-PpoI	PstI	Sse8387I
AscI	BsrGI	MluI	SfiI	StuI
AvrII	Bst1107I	PacI	SgfI	SwaI
BbuI	Bst98I	PaeR7I	SgrAI	XcmI
BclI	Bsu36I	PflMI	SnaBI	XhoI
BglII	ClaI	PinAI	SpeI	

Table 3. Restriction Enzymes That Cut the pAdVAntageTM Vector 6 or More Times.

_

Note: The enzymes listed in boldface type are available from Promega.

 Promega Corporation
 2800
 Woods
 Hollow Road
 Madison, WI
 53711-5399
 USA

 Toll Free in USA 800-356-9526
 Phone 608-274-4330
 Fax 608-277-2516
 www.promega.com

 Part# TB207
 Printed in USA.
 Revised 11/06

V. Related Products

Product	Size	Cat.#
ProFection [®] Mammalian Transfection		
System – Calcium Phosphate	40 transfections	E1200
ProFection [®] Mammalian Transfection		
System – DEAE-Dextran	40 transfections	E1210
Transfectam [®] Reagent for the		
Transfection of Eukaryotic Cells	1mg	E1231
	0.5mg	E1232
TransFast [™] Transfection Reagent	1.2mg	E2431
Tfx™-20 Reagent	4.8mg	E2391
Tfx™-50 Reagent	2.1mg	E1811
Tfx™ Reagents Transfection Trio	5.4mg	E2400
pSI Mammalian Expression Vector	20µg	E1721
pCI Mammalian Expression Vector	20µg	E1731
pCI-neo Mammalian Expression Vector	20µg	E1841
pTARGET TM Mamalian Expression Vector	20 reactions	A1410

VI. References

- 1. Farrell, P.J. *et al.* (1977) Phosphorylation of initiation factor eIF-2 and the control of reticulocyte protein synthesis. *Cell* **11**, 187–200.
- Groskreutz, D. and Schenborn, E. (1994) Increased gene expression in mammalian cell lines using pAdVAntage[™] DNA as a co-transfectant. *Promega Notes* 48, 8–12.
- Kitajewski, J. et al. (1986) Adenovirus VAI RNA antagonizes the antiviral action of interferon by preventing activation of the interferon-induced eIF-2 alpha kinase. Cell 45, 195–200.
- O'Malley, R. et al. (1986) A mechanism for the control of protein synthesis by adenovirus VA RNAI. Cell 44, 391-400.



© 1995-2006 Promega Corporation. All Rights Reserved.

ProFection and Transfectam are registered trademarks of Promega Corporation. pAdVAntage, pTARGET, TransFast and Tfx are trademarks of Promega Corporation.

DNASTAR is a registered trademark of DNASTAR, Inc. GenBank is a registered trademark of US Dept of Health and Human Services.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All prices and specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.

 Promega Corporation
 2800
 Woods
 Hollow Road
 Madison, WI
 53711-5399
 USA

 Toll Free in USA 800-356-9526
 Phone 608-274-4330
 Fax 608-277-2516
 www.promega.com

 Part# TB207
 Printed in USA.
 Revised 11/06