



Promega Points™

Frequent Buyer Program

Collection Card Information

Please Print *(please keep a photocopy of your order until merchandise is received; allow 4-6 weeks for delivery)*

Name: _____

Institution: _____

Department: _____ Room: _____

We cannot deliver to a Post Office Box.

Address: _____

City: _____ State: _____ Zip Code: _____

Phone: _____ Fax: _____

E-mail Address: _____

Ship to Address if different than above:

We cannot deliver to a Post Office Box.

Privacy Notice: Supplying details above gives Promega permission to contact you at each address provided.

Mail Completed Card
With Redemption Cards To:

**PROMEGA CORPORATION
DIRECT MAIL
2800 WOODS HOLLOW RD.
MADISON, WI 53711-5399**

Part #	Item	Pts.	
APPAREL			
A.	PROP155 PH.D. T-shirt (S)	30	<input type="checkbox"/>
	PROP155A PH.D. T-shirt (M)	30	<input type="checkbox"/>
	PROP155B PH.D. T-shirt (L)	30	<input type="checkbox"/>
	PROP155C PH.D. T-shirt (XL)	30	<input type="checkbox"/>
	PROP155D PH.D. T-shirt (XXL)	30	<input type="checkbox"/>
B.	PROP130A Men's Microfleece, 1/4 Zip (S)	80	<input type="checkbox"/>
	PROP130B Men's Microfleece, 1/4 Zip (M)	80	<input type="checkbox"/>
	PROP130C Men's Microfleece, 1/4 Zip (L)	80	<input type="checkbox"/>
	PROP130D Men's Microfleece, 1/4 Zip (XL)	80	<input type="checkbox"/>
	PROP130E Men's Microfleece, 1/4 Zip (XXL)	80	<input type="checkbox"/>
C.	PROP131A Women's Microfleece, 1/4 Zip (S)	80	<input type="checkbox"/>
	PROP131B Women's Microfleece, 1/4 Zip (M)	80	<input type="checkbox"/>
	PROP131C Women's Microfleece, 1/4 Zip (L)	80	<input type="checkbox"/>
	PROP131D Women's Microfleece, 1/4 Zip (XL)	80	<input type="checkbox"/>
	PROP131E Women's Microfleece, 1/4 Zip (XXL)	80	<input type="checkbox"/>
D.	PROP133A Men's Columbia Jacket (S)	100	<input type="checkbox"/>
	PROP133B Men's Columbia Jacket (M)	100	<input type="checkbox"/>
	PROP133C Men's Columbia Jacket (L)	100	<input type="checkbox"/>
	PROP133D Men's Columbia Jacket (XL)	100	<input type="checkbox"/>
	PROP133E Men's Columbia Jacket (XXL)	100	<input type="checkbox"/>
E.	PROP115A Women's Columbia Jacket (S)	100	<input type="checkbox"/>
	PROP115B Women's Columbia Jacket (M)	100	<input type="checkbox"/>
	PROP115C Women's Columbia Jacket (L)	100	<input type="checkbox"/>
	PROP115D Women's Columbia Jacket (XL)	100	<input type="checkbox"/>

LAB EQUIPMENT

F.	PROP095 Waterbath Float	5	<input type="checkbox"/>
G.	PROP053 Fluorescent Ruler	10	<input type="checkbox"/>
H.	A1071 Eluator™ Vacuum Elution Device (set of 4)	40	<input type="checkbox"/>
I.	PROP134 Lab Markers, Black (set of 6)	40	<input type="checkbox"/>
J.	PROP128A Lab Coat (S)	50	<input type="checkbox"/>
	PROP128B Lab Coat (M)	50	<input type="checkbox"/>
	PROP128C Lab Coat (L)	50	<input type="checkbox"/>
	PROP128D Lab Coat (XL)	50	<input type="checkbox"/>
	PROP128E Lab Coat (XXL)	50	<input type="checkbox"/>
K.	PROP135 Lab Markers, Multi-color (set of 8)	60	<input type="checkbox"/>
L.	A7231 Vac-Man® Laboratory Vacuum Manifold	75	<input type="checkbox"/>
M.	PROP098 4-Way Digital Timer	75	<input type="checkbox"/>
N.	PROP052 Benchtop Cooler	100	<input type="checkbox"/>
O.	PROP051 Mini Centrifuge	180	<input type="checkbox"/>
P.	PROP050 Desktop Microcentrifuge	525	<input type="checkbox"/>
Q.	Free Product Credit	Please Inquire	

Part #	Item	Pts.	
OFFICE ACCESSORIES			
R.	PROP100 Mood Magnet	5	<input type="checkbox"/>
S.	PROP126 Post-it® Note Cube (450 sheets)	10	<input type="checkbox"/>
T.	PROP001 Squiggle Clip Pen (set of 3)	10	<input type="checkbox"/>
U.	PROP103 Stress Ball	10	<input type="checkbox"/>
V.	PROP153 Survival Tool	15	<input type="checkbox"/>
W.	PROP073 Desktop Calculator	20	<input type="checkbox"/>
X.	PROP113 Flash Drive (1G)	40	<input type="checkbox"/>
Y.	PROP055 Tool Kit	40	<input type="checkbox"/>
Z.	PROP057 Wireless Presentation/Laser Pointer	60	<input type="checkbox"/>
AA.	PROP099 AM/FM/CD Boombox	100	<input type="checkbox"/>

TRAVEL AND RECREATION

BB.	PROP154 Thermo Tote	20	<input type="checkbox"/>
CC.	PROP104 Water Bottle (20 oz)	20	<input type="checkbox"/>
DD.	PROP006 Soft Pack Insulated Cooler	20	<input type="checkbox"/>
EE.	PROP021 Umbrella	30	<input type="checkbox"/>
FF.	PROP108 Insulated Travel Mug (16 oz)	30	<input type="checkbox"/>
GG.	GE100A Two Movie Tickets* w/Refreshments	30	<input type="checkbox"/>
	<small>* Visit www.hollywoodmoviemoney.com to find a participating theater near you.</small>		
HH.	GE100B Four Movie Tickets* w/Refreshments	45	<input type="checkbox"/>
	<small>* Visit www.hollywoodmoviemoney.com to find a participating theater near you.</small>		
II.	PROP022 Canvas Duffel Bag	80	<input type="checkbox"/>
JJ.	PROP127 Sling Back Pack	80	<input type="checkbox"/>
KK.	PROP112 iPod Shuffle (4GB, 1,000 songs)	150	<input type="checkbox"/>
LL.	PROP116 Portable DVD Player	180	<input type="checkbox"/>
MM.	PROP129 Digital Camera with Case and Memory Card	180	<input type="checkbox"/>

MAGAZINE SUBSCRIPTIONS (1Year) / GIFT CARDS

NN.	Mental Floss	40	<input type="checkbox"/>
OO.	Time Magazine	60	<input type="checkbox"/>
PP.	iTunes Gift Card (\$25)	60	<input type="checkbox"/>
QQ.	Amazon Gift Card (\$25)	60	<input type="checkbox"/>
RR.	National Geographic	90	<input type="checkbox"/>
SS.	Mrs. Field's 48 Nibblers® Bite-size Asst. Cookies	90	<input type="checkbox"/>
TT.	Nature	250	<input type="checkbox"/>
UU.	Science	250	<input type="checkbox"/>

• Please allow 4–6 weeks for delivery of merchandise.

- Items available may change dependent on vendor inventory. Please visit our Web site at: www.promega.com/us/ppoints for the most current list of items available. Promega reserves the right to substitute a similar item if your choice is not available. All photographs are representative samples of actual items. Color and sizes may vary.
- Void where prohibited by law or by institutional policy. Where prohibited from receiving a personal item, a selection of lab equipment and journal subscriptions are available.
- Promega reserves the right to discontinue the program at any time without prior notice.
- We recommend you keep a photocopy of your order until merchandise is received.



Promega