

# Lifesensors, Inc.

Patent Report by Invention

Printed: 11/25/2019 Page 1

COUNTRY	REFERENCE#	TYPE	FILED	SERIAL#	ISSUED	PATENT#	STATUS
---------	------------	------	-------	---------	--------	---------	--------

## BIOLUMINESCENT DETECTION OF PROTEASE ACTIVITY

Inventors: Orcutt, Steven; Yan, Yilin; Wu, Jian; Strickler, James; Butt, Tauseff R.

UNITED STATES	P05084US00	NEW	6/11/2010	61/353,956			EXPIRED
UNITED STATES	P05084US01	FCA	6/10/2011	13/157,734			ABANDONED

## COMPOSITIONS AND METHODS FOR GENDER SORTING

Inventors: Butt, Tauseff R.; Tran, Hiep

UNITED STATES	P02548US0	NEW	4/23/2001	60/286,010			EXPIRED
AUSTRALIA	P02548AU0	DCA	4/23/2002	2002258903			ABANDONED
CANADA	P02548CA0	DCA	4/23/2002	2,445,162	4/5/2011	2,445,162	ABANDONED
EUROPEAN PATENT	P02548EP0	DCA	4/23/2002	02728881.0			ABANDONED
WIPO	P02548WO0	CEQ	4/23/2002	PCT/US02/12590			NAT PHASE
UNITED STATES	P02548US1	DCA	4/23/2004	10/475,433	8/1/2006	7,083,941	ABANDONED

## COMPOSITIONS, METHODS AND KITS FOR ENHANCING PROTEIN EXPRESSION, SOLUBILITY AND ISOLATION

Inventors: Butt, Tauseff R.; Sterner, Savid; Zuo, Xun

UNITED STATES	P04803US00	NEW	12/30/2004	60/640,107			EXPIRED
EUROPEAN PATENT	P04803EP00	DCA	12/30/2005	05855723.2			ABANDONED
WIPO	P04803WO00	CEQ	12/30/2005	PCT/US05/47210			NAT PHASE
UNITED STATES	P04803US01	DCA	5/18/2009	11/794,532	2/21/2012	8,119,369	ISSUED

## DI- AND POLY-UBIQUITIN DEUBIQUITINASE SUBSTRATES AND USES THEREOF

Inventors: Chelur, Dattananda; Eddins, Michael; Leach, Craig; Orcutt, Steven; Peroutka, Raymond J.; Strickler, James; Yan, Yilin

UNITED STATES	P05097US00	NEW	6/24/2010	61/358,177			EXPIRED
UNITED STATES	P05097US01	FCA	6/24/2011	13/168,073	8/27/2013	8,518,660	ISSUED
UNITED STATES	P05097US02	CON	8/26/2013	13/975,941	7/5/2016	9,382,574	ISSUED

## HORMONE RESPONSE SYSTEM

Inventors: Walfish, Paul

UNITED STATES	P02365US0	NEW	8/15/1997	60/055,797			EXPIRED
UNITED STATES	P02365US1	FCA	8/14/1998	09/134,651			ABANDONED
UNITED STATES	P02365US2	CON	9/29/2000	09/677,005			TRANSFER

## METHODS AND COMPOSITIONS FOR ENHANCED PROTEIN EXPRESSION AND PURIFICATION

Inventors: Butt, Tauseff R.; Panavas, Tadas; Karwa, Amolkumar; Peroutka, Raymond J.; Marblestone, Jeffrey G.

UNITED STATES	P04249US00	NEW	12/29/2006	60/877,914			EXPIRED
AUSTRALIA	P04249AU00	DCA	12/28/2007	2007342011			ABANDONED
CANADA	P04249CA00	DCA	12/28/2007	2,674,304	2/23/2016	2,674,304	ISSUED
SWITZERLAND	P04249CH00	DCA	12/28/2007	07866074.3	3/11/2015	2,109,673	ISSUED
CHINA	P04249CN00	DCA	12/28/2007	200780051492.5	4/9/2014	ZL200780051492.5	ISSUED
GERMANY	P04249DE00	DCA	12/28/2007	07866074.3	3/11/2015	60 2007 040 616.1	ISSUED
EUROPEAN PATENT	P04249EP00	DCA	12/28/2007	07866074.3	3/11/2015	2,109,673	NAT PHASE
FRANCE	P04249FR00	DCA	12/28/2007	07866074.3	3/11/2015	2,109,673	ISSUED
UNITED KINGDOM	P04249GB00	DCA	12/28/2007	07866074.3	3/11/2015	2,109,673	ISSUED
INDIA	P04249IN00	DCA	12/28/2007	4399/DELNP/2009	6/29/2017	284769	ISSUED

COUNTRY	REFERENCE#	TYPE	FILED	SERIAL#	ISSUED	PATENT#	STATUS
<i>METHODS AND COMPOSITIONS FOR ENHANCED PROTEIN EXPRESSION AND PURIFICATION continued . . .</i>							
JAPAN	P04249JP00	DCA	12/28/2007	2009-544286	2/7/2014	5469786	ISSUED
SOUTH KOREA	P04249KR00	DCA	12/28/2007	10-2009-7016018	5/29/2015	1526019	ABANDONED
WIPO	P04249WO00	CEQ	12/28/2007	PCT/US07/89035			NAT PHASE
UNITED STATES	P04249US01	FCA	6/1/2010	12/521,468	4/8/2014	8,691,958	ISSUED

### METHODS AND COMPOSITIONS FOR IDENTIFYING LIGANDS FOR NUCLEAR RECEPTORS

Inventors: Tran, Hiep; Askari, Hossein; Schwartz, Michael; Butt, Tauseff R.

UNITED STATES	P02124US0	NEW	2/18/2000	60/183,393			EXPIRED
EUROPEAN PATENT	P02124EP0	DCA	2/20/2001	01912857.8			ABANDONED
WIPO	P02124WO0	CEQ	2/20/2001	PCT/US01/05429			NAT PHASE
UNITED STATES	P02124US1	DCA	12/2/2002	10/204,169	5/2/2006	7,037,656	ABANDONED

### METHODS AND COMPOSITIONS FOR PROTEIN EXPRESSION AND PURIFICATION

Inventors: Butt, Tauseff R.; Weeks, Stephen

UNITED STATES	P02792US0	NEW	1/7/2002	60/346,449			EXPIRED
AUSTRALIA	P02792AU00	DCA	1/7/2003	2003205053	1/29/2009	2003205053	ABANDONED
BELGIUM	P02792BE00	DCA	1/7/2003	03703721.5	12/12/2007	1,470,236	ABANDONED
CANADA	P02792CA00	DCA	1/7/2003	2,472,477	3/20/2012	2,472,477	ISSUED
SWITZERLAND	P02792CH00	DCA	1/7/2003	03703721.5	12/12/2007	1,470,236	ABANDONED
CHINA	P02792-CN	DCA	1/7/2003	N/A			ABANDONED
GERMANY	P02792DE00	DCA	1/7/2003	03703721.5	12/12/2007	603 18 014.0-08	ABANDONED
EUROPEAN PATENT	P02792-EP	DCA	1/7/2003	03703721.5	12/12/2007	1,470,236	ISSUED
FRANCE	P02792FR00	DCA	1/7/2003	03703721.5	12/12/2007	1,470,236	ABANDONED
UNITED KINGDOM	P02792GB00	DCA	1/7/2003	03703721.5	12/12/2007	1,470,236	ABANDONED
IRELAND	P02792IE00	DCA	1/7/2003	03703721.5	12/12/2007	1,470,236	ABANDONED
JAPAN	P02792JP00	DCA	1/7/2003	2003-557532			ABANDONED
UNITED STATES	P02792US1	FCA	1/7/2003	10/338,411	6/13/2006	7,060,461	ISSUED
WIPO	P02792WO0	CEQ	1/7/2003	PCT/US03/00436			NAT PHASE
UNITED STATES	P02792US2	CIP	3/14/2003	10/389,640	5/22/2007	7,220,576	ISSUED
UNITED STATES	P02792US3	DIV	2/28/2006	11/363,660	10/26/2010	7,820,384	ISSUED
UNITED STATES	P02792US4	CON	5/21/2007	11/751,441			ABANDONED

### METHODS AND COMPOSITIONS FOR PROTEIN EXPRESSION AND PURIFICATION

Inventors: Butt, Tauseff R.; Malakhova, Oxana; Malakhov, Michael

UNITED STATES	P03237US00	NEW	6/26/2003	60/482,817			EXPIRED
AUSTRALIA	P03237AU00	DCA	6/28/2004	2004254600			ABANDONED
CANADA	P03237CA00	DCA	6/28/2004	2,530,530			ABANDONED
EUROPEAN PATENT	P03237EP0	DCA	6/28/2004	04756292.1			ABANDONED
INDIA	P03237IN00	DCA	6/28/2004	6050/DELNP/2005			ABANDONED
JAPAN	P03237JP00	DCA	6/28/2004	2006-517749			ABANDONED
WIPO	P03237WO0	CEQ	6/28/2004	PCT/US04/20778			NAT PHASE
UNITED STATES	P03237US1	DCA	11/15/2005	10/504,785	2/2/2010	7,655,413	ISSUED

### SYNTHETIC PROTEASE SUBSTRATES, ASSAY METHODS USING SUCH SUBSTRATES AND KITS FOR PRACTICING THE ASSAY

Inventors: Butt, Tauseff R.; Manimala, Joseph; Cejas, Mabel; Strickler, James; Kingsbury, William

UNITED STATES	P04879US00	NEW	2/3/2010	61/300,987			EXPIRED
---------------	------------	-----	----------	------------	--	--	---------

COUNTRY	REFERENCE#	TYPE	FILED	SERIAL#	ISSUED	PATENT#	STATUS
<i>SYNTHETIC PROTEASE SUBSTRATES, ASSAY METHODS USING SUCH SUBSTRATES AND KITS FOR PRACTICING</i>							
CANADA	P04879CA00	CEQ	2/3/2011	2,730,618			ABANDONED
UNITED STATES	P04879US01	FCA	2/3/2011	13/020,416	4/15/2014	8,697,376	ISSUED

END OF REPORT

TOTAL ITEMS SELECTED = 64